



DIANTHUS uHTS
PRODUCT MANUAL

What is included in this product manual

DIANTHUS uHTS

User Manual

Table of contents

1.	About this user manual	6	3.2.1.	Technical specifications	11
1.1.	Directions for more detailed information	6	3.2.2.	Connections for input and output	12
2.	Safety information	7	3.2.3.	Replacement of fuses	13
2.1.	Symbols and descriptions	7	3.2.4.	LED panel and device status	13
2.2.	Use and misuse	8	3.2.5.	Stacking of instruments	13
2.3.	Safety instructions	8	3.3.	Legal	14
3.	The Dianthus uHTS system	9	3.4.	Limited warranty	14
3.1.	General	9	4.	Dianthus uHTS setup	15
3.1.1.	Intended use	9	4.1.	Scope of delivery	15
3.1.2.	Conformity	9	4.2.	Unpacking	15
3.1.3.	Identification	10	4.3.	Startup	15
3.2.	Technical information	11	4.4.	Cleaning	16

4.5.	Software updates	16	5.8.	Data storage and export	20
5.	Using Dianthus uHTS	16	6.	Troubleshooting	20
5.1.	Dianthus uHTS well plates	16	6.1.	Wrong insertion of Dianthus uHTS well plates	20
5.1.1.	General tips for handling plates	17	6.2.	Disconnect from control PC	20
5.2.	Barcode Reading	17	6.3.	Restarting Dianthus uHTS	20
5.3.	Dianthus uHTS Control software	17	6.4.	Customer support	20
5.4.	Underlying physical principles	18	7.	Patents and intellectual property	21
5.5.	How does a Dianthus uHTS measurement look like?	18	8.	Transport and disposal	21
5.6.	Sequence of positioning and data acquisition for one well in a 384-microwell plate	19	8.1.	Repackaging for transport	21
5.7.	Sequence of positioning and data acquisition for a 1536-microwell plate	20	8.2.	Waste disposal	21
			8.3.	System disposal	21

1. About this user manual

This user manual gives guidance on the correct use of the Dianthus uHTS system. It covers system specifications, safety considerations and installation as well as why and how to run experiments with Dianthus uHTS. Please read this manual carefully before starting and make sure the contents are fully understood. Keep this manual available near the system for future reference. In case of loss, please visit the NanoTemper Technologies Support Center (<https://support.nanotempertech.com/hc/en-us>) for a replacement copy of this manual.



1.1. Directions for more detailed information

For further, more detailed information on scientific principles and recommendations for assay development as well as software usage visit our Support Center (<https://support.nanotempertech.com/hc/en-us>).

The NanoTemper Technologies **Support Center** is an online community to obtain resources for NanoTemper products such as application-notes, tech-notes or protocols. You can also pose questions and have them answered by NanoTemper support guides

Do not miss out on our online-tool “Degree of labeling calculator” that helps with calculation of the degree-of-labeling (DOL): <https://nanotempertech.com/user-tools/dol-calculator/>

2. Safety information

To ensure operation safety, this system must be operated correctly. Carefully read this chapter to fully understand all necessary safety precautions before operating the system.

2.1. Symbols and descriptions

This section describes the safety symbols and descriptions used in this manual, as well as the labels on the system.

Please take a moment to understand what the signal words **WARNING!** **CAUTION** and **NOTE** mean in this manual.

WARNING! A **WARNING!** indicates a potentially hazardous situation which, if not avoided, could result in serious injury or even death.

CAUTION A **CAUTION** indicates a potentially hazardous situation which, if not avoided, may result in minor or moderate injury. **CAUTION** may also be used to alert against damaging the equipment or the system.

Do not proceed beyond a **WARNING!** or **CAUTION** notice until you understand the hazardous conditions and have taken the appropriate steps.

NOTE A **NOTE** provides additional information to help the operator achieve optimal system and assay performance.



Read manual label. This label indicates that you must read the manual before using the system. This label is positioned on the back of the instrument.



Warning symbol. This symbol, when used on its own or in conjunction with any of the following icons indicates the need to consult the provided manual, because a potential risk exists if the operating instructions are not followed. This label is positioned on the back of the instrument.



Warning symbol. This symbol indicates moving parts that can crush and cut. This warning label is positioned on the instrument tray holding the microwell plate.



Warning symbol. This symbol indicates the presence of electric shock hazards. To avoid risk of injury from electric shock, do not open the enclosure. The enclosure should only be opened by NanoTemper authorized personnel. This label is positioned on the back of the device.

2.2. Use and misuse

Use the Dianthus uHTS system only after having read and fully understood this user manual. Use the system only in perfect condition. If the system shows any signs of damage, stop operation and contact NanoTemper Technologies customer support.

Do not modify the system in any way. Do not use it for anything other than its intended purpose.

2.3. Safety instructions

WARNING! The door of the system can pinch or injure your hands or fingers. Keep fingers safe while opening and closing the door. Do not touch the instrument while parts are moving. Do not reach into the opening when the door is open.

WARNING! Connect the Dianthus uHTS to the AC power supply using the supplied power cable. Since the instrument is assembled in line with the specifications for safety class IEC 61010-1:2010, it must only be connected to an outlet that has a ground contact.

WARNING! Danger of electric shock, fire and skin burns. Do not open the system (other than operating the door/loading hatch via the software). Do not reach into the door opening.

WARNING! Using hazardous or infectious substances in the system may pose a risk of explosion, implosion, release of gases or infection. Use only non-hazardous, non-infectious, aqueous samples. Dispose of used microwell plates according to the substances contained in them and according to locally applicable regulations concerning chemical waste.

CAUTION The instrument contains an IR-laser module (invisible laser radiation class 3B according to IEC 60825-1: 2014) and a green laser module (visible laser radiation class 3B according to IEC 60825-1: 2014). Lasers or laser systems emit intense, coherent electromagnetic radiation that has the potential of causing irreparable damage to human skin and eyes. Direct eye contact can cause corneal burns, retinal burns, or both, and possible blindness. Do not attempt to open the instrument as this poses a risk of personal injury or damage to the instrument.

When the instrument is used as intended it protects the user from dangerous laser radiation and is classified as a **class 1 Laser Product**.



CAUTION The system must be installed in a way that does not hinder access to the power switch and power plug.

CAUTION Do not replace the detachable power cord with an inadequately rated cord.

CAUTION Broken glass can cut skin. Do not use if the front or back glass is broken.

CAUTION The weight of the Dianthus uHTS instrument is approx. 70 kg, do not move the instrument alone (four persons required for transport). If you move the instrument alone, you risk personal injury or damage to the instrument.

CAUTION Only NanoTemper Technologies staff may service and open the instrument.

CAUTION Disconnect the mains supply before changing the fuses.

CAUTION This equipment is not intended for use in residential environments and may not provide adequate protection to radio reception in such environments. According to IEC 61326-1, in connection with CISPR 11, this device is: Group 1, Class A

Group 1: Equipment has intentionally generated and/or uses conductively-coupled radio frequency energy that is necessary for the internal function of the equipment itself.

Class A: Equipment is suitable for use in all establishments other than domestic and those directly connected to a low-voltage power supply network that supplies buildings used for domestic purposes. There may be potential difficulties in ensuring electromagnetic compatibility in other environments due to conducted and radiated disturbances.

NOTE Insufficient air supply can cause overheating of the system. Assure enough air supply by not covering the back of the system. Leave at least 15 cm of space between system and any wall or other obstruction.

3. The Dianthus uHTS system

3.1. General

3.1.1. Intended use

The Dianthus uHTS system provides fast and highly sensitive detection and quantification of molecular interactions in microwell plates. The system is intended for research purposes only. It is not to be used for diagnostic purposes.

3.1.2. Conformity

The following safety and electromagnetic standards were considered:

- IEC 61010-1:2010/AMD1:2016, Safety requirements for electrical equipment for measurement, control and laboratory use.
- IEC 61326-1:2021 EMC, Electrical equipment for measurement, control and laboratory use – EMC requirements.
- IEC 60825-1:2014, Safety of laser products.
- Complies with 21 CFR 1040.10 and 1040.11 except for conformance with IEC 60825-1 Ed. 3., as described in Laser Notice No. 56, dated May 8, 2019

3.1.3. Identification

The identification label (Figure 1) is positioned at the rear panel of the instrument. It includes manufacturer information, system model name and serial number (SN), electrical requirements, and the CE conformity symbol.

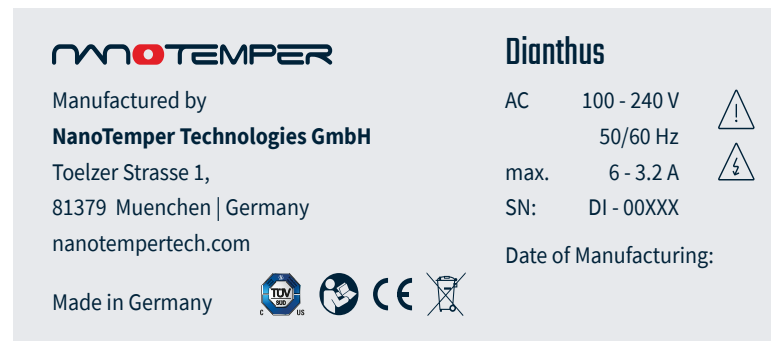


Figure 1: Identification label for Dianthus uHTS.

3.2. Technical information

3.2.1. Technical specifications

Electricity

Input Voltage AC	AC 100-240 V -10 % +10 %
Voltage phase	Single phase
Mains frequency	50/60 Hz
Overvoltage category	CAT II
Input current AC	6 - 3.2 A
Pollution degree	2
Fuse	Fuse link 5 x 20 mm, 10 A, 250 V, time-lag T (2x)

Environmental

Operating temperature	20 – 30 °C (indoor use only)
Storage temperature	-20 – 30 °C
Humidity	5–70 %, non-condensing
Operating altitude	Max. 2000 m

Dianthus uHTS dimensions

Width	61 cm (24.0")
Height	42 cm (16.5")

Depth

57 cm (22.4"), with open tray: 69 cm (27.2")

Weight

70 kg (154.3 lbs) net

Green Laser

Wavelength	520 nm ± 10 nm
Power	10 mW max.

IR Laser

Wavelength	1475 nm ± 15 nm
Power	120 mW max.

Dianthus uHTS Laser classification

The device is LASER PRODUCT CLASS 1

Temperature control

Temperature control range	20 °C – 25 °C
Maximum difference to room temperature*	± 5 °C
Precision of temperature control	± 0.25 °C

*for best possible results we recommend operating Dianthus uHTS at a temperature setpoint equal to the ambient temperature of the room.

3.2.2. Connections for input and output

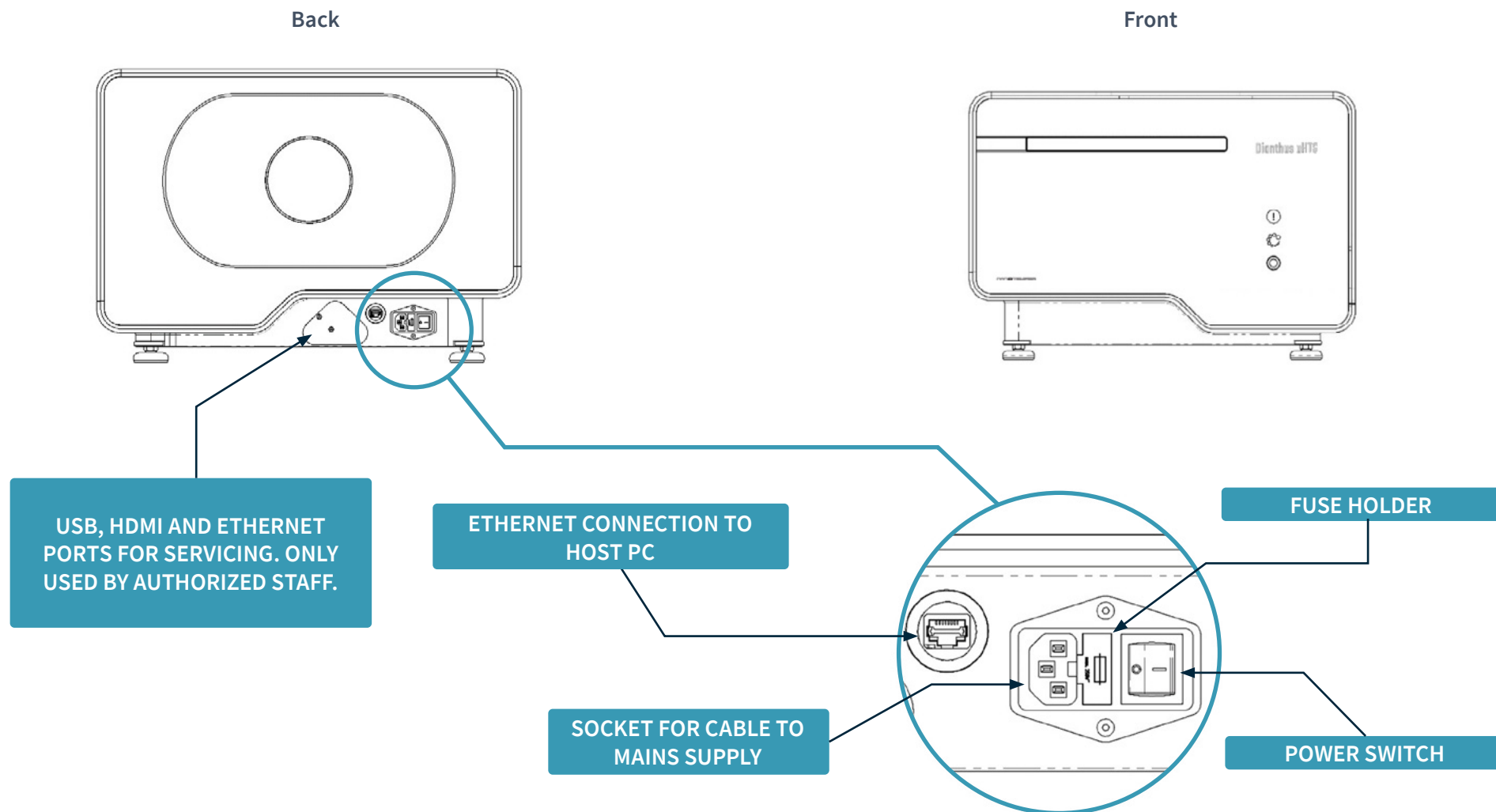


Figure 2: Connections on the Dianthus uHTS device.

Type	Function	Position
Ethernet	To connect the system to the Control PC/LAN via Ethernet cable.	Back panel
AC Power	To connect the system to electrical power.	Back panel

All connections can be found at the rear of the instrument.

3.2.3. Replacement of fuses

We recommend not to change the fuses on your own. Please consult NanoTemper Support (<https://nanotempertech.com/support>) if you think the fuse is broken. To replace the fuses, disconnect the mains supply and open the fuse holder (see Figure 2). Replace the two fuses with new ones of type fuse link 5 x 20 mm, 10 A, 250 V, time-lag T (available, e.g. from ESKA, art. no. 522727).

3.2.4. LED panel and device status



Warning state: Appears when there is an error state, such as a failed initialization. Open DI.Control to receive more detailed information about the device status. Contact support if issue cannot be resolved by power-cycling the instrument.



Measuring state: Indicates that the instrument is actively measuring and acquiring data. Even if connection to the control PC is lost or interrupted, measurement continues, and data is saved.



Initialization: Flashes yellow while initializing and while temperature setpoint is not yet reached. Blue when instrument is initialized and ready.

3.2.5. Stacking of instruments

To save space in your lab, it is possible to stack a maximum number of two instruments on top of each other.

3.3. Legal

1. NanoTemper Technologies shall not be held liable, either directly or indirectly, for any consequential damage incurred as a result of product use.
2. Prohibitions on the use of NanoTemper Technologies software:
 - Copying software for purposes other than backup
 - Transfer or licensing of the right to use software to a third party
 - Disclosure of confidential information regarding software
 - Modification of software
 - Use of software on multiple workstations, network terminals, or by other methods
3. The content of this manual is subject to change without notice for product improvement.
4. This manual is considered complete and accurate at publication.
5. This manual does not guarantee the validity of any patent rights or other rights.
6. If a NanoTemper Technologies software program has failed, causing an error or improper operation, this may be caused by a conflict from another program operating on the controlling PC. In this case, take corrective action by uninstalling the conflicting product(s).
7. NanoTemper and Dianthus are registered trademarks of NanoTemper Technologies GmbH in the United States of America and other countries.
8. Unauthorized resale is not permitted.

3.4. Limited warranty

Products sold by NanoTemper Technologies, unless otherwise specified, are warrantied to be free of defects in materials and workmanship for a period of one year from the date of shipment. If any defects in the product are identified during this warranty period, NanoTemper Technologies will repair or replace the defective part(s) or product free of charge.

This warranty does not apply to defects resulting from the following:

1. Improper or inadequate installation.
2. Improper or inadequate operation, maintenance, adjustment or calibration.
3. Unauthorized modification or misuse.
4. Use of unauthorized microwell plates and accessories.
5. Use of consumables, disposables and parts not supplied by an authorized NanoTemper Technologies distributor.
6. Corrosion due to the use of improper solvents, samples, or due to surrounding gases.
7. Accidents beyond NanoTemper Technologies' control, including natural disasters.

This warranty does not cover consumables like microwell plates, reagents, labeling kits and the like. It also does not cover normal wear-and-tear.

The warranty for all parts supplied and repairs provided under this warranty expires on the warranty expiration date of the original product. For inquiries concerning repair service, contact NanoTemper Technologies after confirming the model name and serial number of your NanoTemper Technologies system (see 3.1.3).

4. Dianthus uHTS setup

The Dianthus uHTS should be installed by NanoTemper Technologies personnel to ensure safety measures are taken and to confirm proper functionality of the instrument.

4.1. Scope of delivery

Upon receiving the system, please check package contents for completeness. The Dianthus uHTS system package contains the following items:

Item	Description
Dianthus uHTS system	-
User manual	This user manual
Cables	Power cord for power supply, Network cable for connection to control PC
Control PC and Monitor	Desktop control PC with monitor for Dianthus uHTS system

4.2. Unpacking

The Dianthus uHTS system should only be unpacked and installed by trained NanoTemper Technologies personnel to ensure proper functionality of the instrument upon delivery.

Note: After transport it is possible that the instrument is cool and needs some time to equilibrate to room temperature. If the instrument is turned on while the instrument temperature is significantly lower than the set temperature the measurement accuracy may be reduced.

4.3. Startup

To begin using the Dianthus uHTS system:

- 1. Power Connection:** Connect the Dianthus uHTS system to power by plugging in the power supply cable.
- 2. Connect to Control PC (optional):** Connect the Dianthus uHTS system to the control PC using the ethernet connection located at the back of the instrument.
- 3. Power On:** Switch on the power switch.
- 4. Initialization:** The LED display will indicate the initialization process. Once ready, the LED display will show the “Connected” state, with the initialization LED glowing blue.

Communication with the Dianthus uHTS instrument

The instrument features a Google Remote Procedure Call (gRPC) interface accessible via port number 172.16.4.1:5002. Communication with the instrument’s gRPC interface can be established through two methods, depending on whether the instrument is used in a manual mode or in an automated mode:

- 1. Using the Pre-installed DI.uHTS Control Interface (manual mode):** A control interface, DI.uHTS Control, is pre-installed on the instrument for convenient and straightforward access. We recommend utilizing the Microsoft Edge browser pre-installed on the NanoTemper control PC to access this control interface. Upon launching the browser, the bookmark to the control interface is already set. Note: If you encounter the message “your connection isn’t private”, click “advanced” to grant permissions for uninterrupted access.
- 2. Developing Your Own Driver for gRPC Interface (automated mode):** For full automation and customization, users have the option to develop their own driver for the instrument’s gRPC interface. By creating a custom driver, users can tailor the instrument’s functionalities to suit their specific requirements and seamlessly integrate it into their existing automation workflows.

4.4. Cleaning

We recommend regular maintenance of the instrument by trained NanoTemper Technologies personnel to ensure internal components are functional. To clean the outside surface of the system, unplug the power supply at the back. Wipe the surface, including the front display, with a cloth slightly dampened with water or ethanol.

4.5. Software updates

Embedded System Updates

Software updates for the embedded system are exclusively conducted by authorized NanoTemper Technologies personnel during regular maintenance visits. This ensures that the embedded system remains up-to-date and operates optimally.

DI.uHTS Control Software Updates

Users have the capability to perform software updates for the DI.uHTS Control software independently. This feature empowers users to keep the control software current, ensuring compatibility with evolving requirements and enhancing the overall functionality of the system.

5. Using Dianthus uHTS

The Dianthus uHTS instruments utilize a unique photophysical property of fluorophores to detect and quantify binding events rapidly and effortlessly. This characteristic revolves around a phenomenon known as Spectral Shift, which involves observing changes in the fluorescence emission spectrum. Specifically, Spectral Shift relies on detecting alterations in the emission of a fluorescent probe, employing a ratiometric dual-wavelength method. This approach enables the characterization of interactions and the derivation of affinity constants and is further described in section 5.3.

Measurements in Dianthus uHTS instruments occur in industry standard microwell plates.

5.1. Dianthus uHTS well plates

NOTE Dianthus uHTS instruments can only be used with Dianthus uHTS 384- or 1536-well plates which can be purchased from NanoTemper Technologies.

CAUTION Using any other plate with the instrument could result in damage or unreliable results.

The well geometry and the specific coating of Dianthus uHTS 384- and 1536-well plates ensure highest reproducibility and data quality. The optimal working volume for 384-well plates is 20 μL , while the optimal working volume for 1536-well plates is 7 μL . Both plate types are coated with a special polymer to avoid protein adsorption to well walls and well bottom. The bottom of the plate is made of transparent material as the Dianthus uHTS optical system measures the sample from the bottom. This has the added advantage that the top of the plate can be sealed to avoid sample oxidation and evaporation.

5.1.1. General tips for handling plates

- Do not leave fingerprints on the foil bottom. In case, wiping the bottom with a lint-free tissue can remove fingerprints.
- Avoid dust and scratches on the plate bottom. Importantly, one should avoid taking too much dust into the instrument. In case, wiping the bottom with a lint-free tissue can remove dust.
- Use reverse pipetting where possible to avoid air bubbles. If you pipette manually in a 384-well plate it is advisable to use reverse pipetting in cases where it is possible (e.g. when dispensing the buffer for a dilution series).
- In general, avoid air bubbles at the plate bottom. This can affect the measurement. If air bubbles appear, centrifuge the plate **open** (i.e. without seal).
- For few samples that are pipetted manually, first pipette in tubes and then transfer to the plate using reverse pipetting. In tubes it is easier to see while pipetting, therefore results are often more accurate.
- When pipetting directly and manually in the plate, avoid piercing the foil bottom with your pipette tip.
- One should always work with the optimal working volume of 20 μ L per well for 384-well plates and 7 μ L per well for 1536-well plates.
- If plates are kept in a fridge or in an incubator with significantly different temperature than the device temperature, it is advisable to incubate the plates at least 30 min before measuring at device temperature. At best, pre-incubation is done inside the instrument.
- Ensuring thorough mixing of the sample prior to measurement is essential for achieving consistent results and maintaining high data quality in Dianthus uHTS measurements.
- For 384-well plates, we advise conducting a minimum of 15 mixing cycles, with the sample occupying at least 80% of the final volume when using liquid

handling solutions with pipette tips. Alternatively, when employing contact-less liquid handling, we recommend shaking the plate at a minimum of 10 G for 180 seconds, followed by a 30-minute incubation period.

- When dispensing nanoliter volumes into 1536-well plates, adding the target solution is typically sufficient for adequate mixing. We recommend adjusting incubation times based on signal stability testing for optimal results.

5.2. Barcode Reading

Microwell plates can be purchased from NanoTemper Technologies with customer-defined barcodes and positions. We support all common barcode formats, e.g. Code 128 or Interleaved 2 of 5 symbology. Plates can be customized with barcode labels at all four edges of the microwell plates, however, one barcode at the right/EAST side is compulsory since the system will only read the barcode on this side.

5.3. Dianthus uHTS Control software

The Dianthus uHTS instrument is accompanied by the DI.uHTS Control software, providing comprehensive control and monitoring capabilities for the instrument's operations.

Key Features:

1. **Instrument Control:** Users can efficiently manage the instrument's settings through the DI.uHTS Control software. This includes selecting the well plate type, specifying the wells to be measured, and adjusting parameters such as LED power and device temperature.
2. **Real-time Feedback:** During measurement, the software offers real-time feedback for each well. This includes indicators for fluorescence signal sufficiency, detector saturation, pre-scan status, well-scan analysis and whether the ratio value surpasses or falls below the predefined threshold.

3. **Ratio Value Visualization:** The software facilitates easy visualization of ratio values for each well. Users can hover over individual wells to access corresponding ratio values, enhancing data interpretation and analysis.
4. **Data Export:** Users have the option to export data from the DI.uHTS Control software in JSON file format. This allows for seamless integration with various analysis tools and platforms, facilitating in-depth data analysis and interpretation.
5. **Data Management:** DI.uHTS Control software allows users to review past experiments through a comprehensive data experiment history. Additionally, users can conveniently export corresponding raw data files for further analysis or archival purposes.

5.4. Underlying physical principles

Spectral Shift detection is a biophysical technique that measures the strength of the interaction between two molecules by detecting a variation in the fluorescence emission profile of a fluorescently labeled or intrinsically fluorescent target when a binding event occurs.

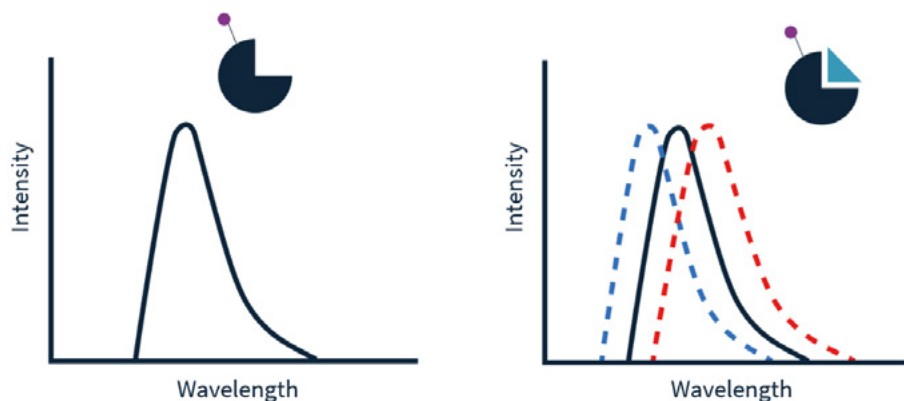


Figure 3: Illustration of a binding-induced change in the fluorescence emission spectrum. Spectrum can shift either to longer wavelengths (red-shift), or to shorter wavelengths (blue-shift).

The emission wavelength is strongly related to the chemical environment of the fluorophore, which can be changed by the binding of a ligand to the target. Chemical environment changes for the fluorophore through ligand binding can occur by 2 distinct mechanisms: either by binding of a ligand in close proximity to the position of the fluorophore or by inducing a conformational change in the target molecule (Figure 4).

5.5. How does a Dianthus uHTS measurement look like?

The Dianthus uHTS instrument accommodates either 384- or 1536-microwell plates as sample vessels, providing flexibility to suit diverse experimental needs. For data acquisition, the instrument applies SpS measurements. A full 384-well plate can be measured in approx. 30 minutes, while a full 1536-well plate requires approx. 7 minutes for completion.

Users have the option to utilize the instrument as a standalone device or seamlessly integrate it into laboratory automation setups using Google Remote Procedure Call (gRPC) interface. This integration capability enables enhanced efficiency and workflow automation. For comprehensive information on automation options, please refer to Section 4.3 of this user manual. Alternatively, feel free to reach out to our support team for personalized assistance.

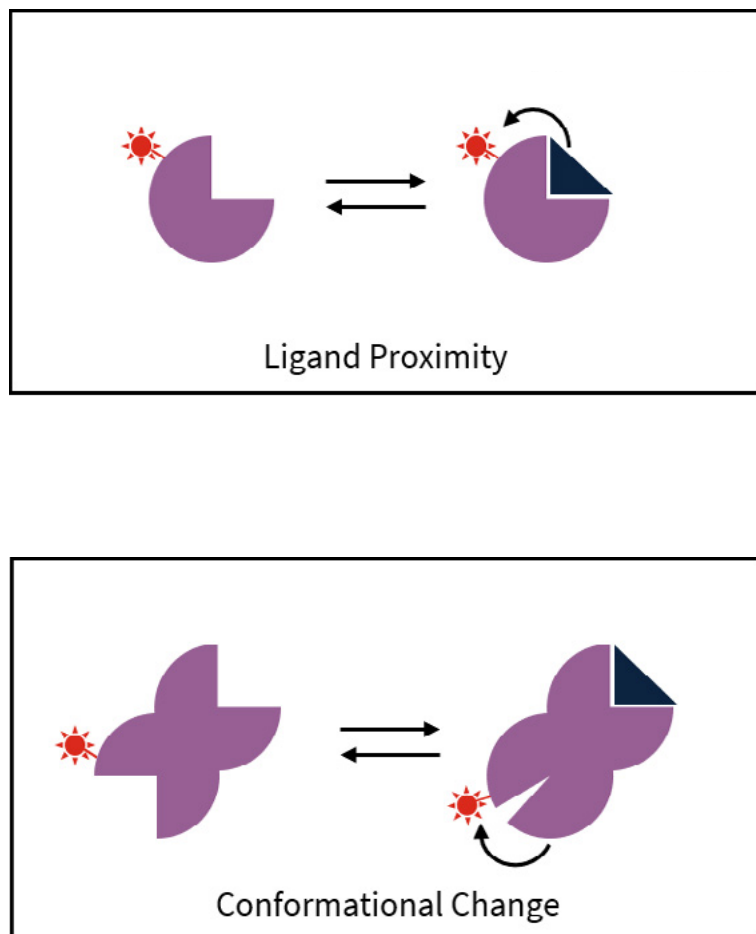


Figure 4: Illustration of chemical environment changes for fluorescent probes upon target-ligand complex formation.

5.6. Sequence of positioning and data acquisition for one well in a 384-microwell plate

Before one well can be measured, the optical system needs to be carefully positioned.

As a first step of optics positioning, Dianthus moves the optical setup in Z-direction and uses reflected laser light to determine the accurate position of the interface

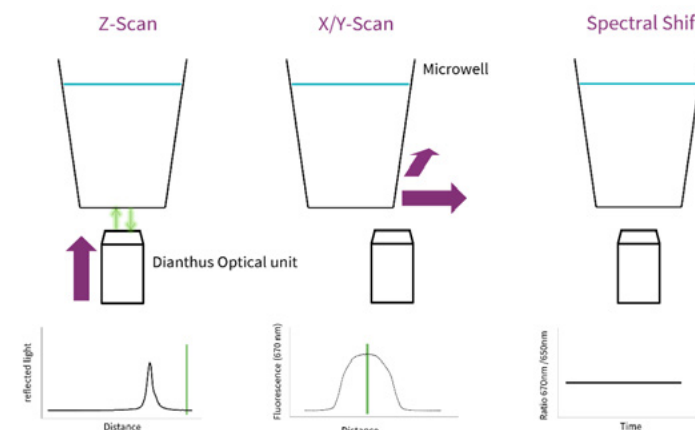


Figure 5: Illustration of the serial steps of establishing the spatial dimensions of one specific well. Z-Scan, X-Scan and Y-Scan and ensure high precision positioning of the optical system before a spectral shift measurement is recorded for that well.

between air and plate bottom, as well as plate-bottom and sample (Figure 5 left). This careful and highly precise positioning of the optical system is pivotal. The Z-scan is followed by a sample scan to measure sample fluorescence while the well moves in X- and Y-direction above the optics (Figure 5 center). From the resulting well profile, Dianthus determines the well center in X and Y, here indicated as a green line. Following optics positioning, Dianthus now performs the spectral shift measurement (Figure 5 right). The Z-position for a spectral shift measurement lies at a distance of 1.5 mm from the well-bottom.

5.7. Sequence of positioning and data acquisition for a 1536-microwell plate

When measuring 1536-microwell plates, the system first carries out a calibration measurement (pre-scan) to determine the accurate X-, Y- and Z-positions of all wells. For this purpose, the above-described Z-scans and X/Y-scans are carried out at selected wells only. To enable the system to measure these positions, it is important that all selected wells contain a fluorescent sample – empty wells cannot be detected. Subsequently, a fast uHTS scan of the entire plate is performed. The entire measurement cycle, including opening the drawer, closing the drawer, calibration and uHTS-scan takes approximately 7 minutes.

5.8. Data storage and export

All measured data are stored on the device and can be accessed and downloaded via the history in the DI.uHTS Control software. When using the DI.uHTS Control software, the user will receive a warning when the memory use reaches 80%. When using the system via a custom driver, the user should query the memory state at regular intervals. When approaching the memory limit, we recommend deleting the oldest data to free memory. The system will never delete data automatically. If the memory limit is reached, an error will be displayed, and all further measurements are blocked.

Generated data can be exported from the DI.uHTS Control software in JSON file format. This allows for seamless integration with various analysis tools and platforms, facilitating in-depth data analysis and interpretation.

For comprehensive guidance on integrating data from the DI.uHTS Control software into external analysis platforms such as Gene Data Screener or Spotfire, we encourage you to contact our support team. Our experts will provide detailed information and assistance to ensure seamless integration and optimal utilization of your data within your preferred analysis environment.

6. Troubleshooting

6.1. Wrong insertion of Dianthus uHTS well plates

In case a plate is inserted in a wrong way and blocks the instrument, the drawer cannot close, the instrument will stop and go into a warning state. In such a case please switch off the instrument and contact the NanoTemper Technologies support (www.nanotempertech.com/support).

6.2. Disconnect from control PC

In case of a disconnect between instrument and control PC the instrument would still complete the ongoing measurement. This is indicated by the LED panel. Please do not switch off the instrument in this state. Wait until the measurement is complete and reestablish the connection. The DI.Control software will then obtain the measured data from the internal instrument PC.

6.3. Restarting Dianthus uHTS

In case the system freezes, wait one minute. If it does not un-freeze, use the switch at the back of the system to switch off the instrument. Wait 30 seconds for complete shutdown, then restart the instrument. The system will start up again automatically.

6.4. Customer support

In case of any issues not described in this user manual, please don't hesitate to contact NanoTemper Technologies customer support at: www.nanotempertech.com/support.

7. Patents and intellectual property

Dianthus™ and NanoTemper® are registered trademarks of NanoTemper Technologies. Products are protected by US8431903B2, US8853650B2, US9459211B2, US10345312B2, US8741570B2 and further patent applications owned by NanoTemper Technologies. Purchase of products related to this technology from NanoTemper Technologies provide an implied license for non-profit and in-house research use. Any application of above-mentioned technologies, e.g. for in-vitro diagnostic and medical use, require a separate license from NanoTemper Technologies. A license may be granted by NanoTemper Technologies on a case-by-case basis and is entirely at NanoTemper Technologies' discretion.

8. Transport and disposal

8.1. Repackaging for transport

The Dianthus uHTS instrument should be repacked only by trained NanoTemper Technologies personnel to ensure safety and stability during transport. Please store the instrument box for that purpose. If the instrument box was discarded NanoTemper Technologies can provide replacement at the cost of packaging material and shipment.

8.2. Waste disposal

Please dispose of used microwell plates according to the substances contained in them and according to locally applicable regulations concerning chemical and glass waste.

8.3. System disposal

The system may need to be decontaminated before disposal. Please contact NanoTemper Technologies for more information.



This symbol indicates that this system may not be disposed of as unsorted municipal waste and must be collected separately. It must be disposed of according to locally applicable regulations regarding electrical and electronic equipment. The symbol is positioned at the back of the instrument.

Contact

NanoTemper Technologies GmbH

Global headquarters:
Toelzer Strasse 1,
81379 Muenchen | Germany

Phone: +49 (0)89 4522895 0

Fax: +49 (0)89 4522895 60

info@nanotempertech.com
nanotempertech.com

Dianthus™ is a trademark, registered with the U.S. Federal Trademark registration.

NanoTemper® is a registered trademark and registered in the U.S. Patent and Trademark Office.

V01_2024-09-06



NanoTemper cares for the future of our planet, therefore we want to do our best to protect it and reduce the environmental footprint of our products.

*This product manual is printed in Germany, using **EU Ecolabel approved paper from certified FSC® responsible forestry.***

A decorative graphic consisting of several thick, blue, irregular lines that flow across the page, creating a modern, organic frame around the text.

NanoTemper Technologies GmbH

Global headquarters:

Toelzer Strasse 1,

81379 Muenchen | Germany

Phone: +49 (0) 89 4522895 0

Fax: +49 (0) 89 4522895 60

info@nanotempertech.com

nanotempertech.com

NanoTemper[®], and Dianthus[™] are registered trademarks.