

## Monolith Protocol M0-P-038

# IDO1 – NLG919 Analogue

IDO1 (Indoleamine 2,3-dioxygenase 1) is a monomeric heme-containing dioxygenase that catalyzes the oxidative cleavage of L-Tryptophan. IDO1 shows elevated expression in some human cancers and is thought to play a role in the tumor immuno-editing process which sets up peripheral tolerance to tumor antigens. It has therefore gained interest as a drug target in cancer therapy, but no inhibitor has advanced to a drug so far. The interaction described in this protocol has been used as a positive control to validate fragments from a screening approach to identify new leads for IDO1 inhibitors.

protein – small molecule interaction | fragment screening

### A1. Target/Fluorescent Molecule

Indoleamine 2,3-dioxygenase 1

[uniprot.org/uniprot/PI4902](https://uniprot.org/uniprot/PI4902)

### A2. Molecule Class/Organism

Dioxygenase

*Homo sapiens (Human)*

### A3. Sequence/Formula

MAHAMENSWT ISKEYHIDEE VGFALPNPQE NLPDFYNDWM FIAKHLPLDI ESGQLRERVE KLNMLSIDHL TDHKSQRLAR  
LVLGCITMAY VWGKGHGDVR KVLPRNIAVP YCQLSKKLEL PPILVYADCV LANWKKKDPN KPLTYENMDV LFSFRDGDSC  
KGFFLVSLLV EIAAASAIKV IPTVFKAMQM QERDTLLKAL LEIASCLEKA LQVFHQIHDH VNPKAFFSVL RIYLSGWKGN  
PQLSDGLVYE GFWEDPKEFA GGSAGQSSVF QCFDVLLGIQ QTAGGGHAAQ FLQDMRRYMP PAHRNFLCSL ESNPSVREFV  
LSKGDAGLRE AYDACVKALV SLRSYHLQIV TKYILIPASQ QPKENKTSER PSKLEAKGTG GTDLMNFKLT VRSTTEKSL  
KEG

### A4. Purification Strategy/Source

Enzyme cloned from human cDNA, expressed in E. coli

Giotto Biotech (Giotto Biotech S.r.l., Firenze)

### A5. Stock Concentration/Stock Buffer

1 mg/mL | 20 µM

20 mM Tris-HCl, pH 8.0, 100 mM NaCl

### A6. Molecular Weight/Extinction Coefficient

45.3 kDa

51,380 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

## A7. Dilution Buffer

50 mM Tris-HCl, pH 7.4, 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 2 mM DTT, 0.05% TWEEN® 20

## A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS (MO-L001, NanoTemper Technologies GmbH)

1\* Labeling Buffer NHS | 1\* A-Column | 1\* Dye RED-NHS 2nd Generation (10 µg) | 1\* B-Column

## A9. Labeling Procedure

1. Add 50 µL of Labeling Buffer NHS containing 0.05% TWEEN® 20 to 50 µL of 20 µM IDO1 to obtain 100 µL of a 10 µM solution.
2. Use the A-Column to perform a buffer exchange into Labeling Buffer NHS.
  - a. Invert A-Column to suspend slurry and twist off bottom (twist slightly in both directions).
  - b. Loosen the cap of the column and place it in a 1.5 mL microcentrifuge collection tube.
  - c. Centrifuge at **1500 × g** for **1 min** to remove excess liquid.
  - d. Add 300 µL of Labeling Buffer NHS and centrifuge at **1500 × g** for **1 min** (3x).
  - e. Place 100 µL of the 10 µM IDO1 solution in the center of the resin.
  - f. Place the sample in a **new** microcentrifuge collection tube and centrifuge at **1500 × g** for **2 min**.

The collected flow-through should yield around 100 µL of ~7.5 µM IDO1 (70 – 80% recovery).
3. Add 30 µL of DMSO to 10 µg RED-NHS dye to obtain a ~470 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 8.5 µL of the 470 µM dye solution with 91.5 µL of Labeling Buffer NHS to obtain 100 µL of a 39.8 µM dye solution (5x protein concentration).
5. Mix IDO1 and dye in a 1:1 volume ratio (200 µL final volume, 5.8% final DMSO concentration).
6. Incubate for 30 minutes at room temperature in the dark.
7. In the meantime, remove the top cap of the B-Column and pour off the storage solution. Remove the bottom cap and place with adapter in a 15 mL tube.
8. Fill the column with dilution buffer and allow it to enter the packed resin bed completely by gravity flow. Discard the flow through collected. Repeat this step 3 more times.
9. Add 200 µL of the labeling reaction from step 5 to the center of the column and let sample enter the bed completely.
10. Add 300 µL of dilution buffer after the sample has entered and discard the flow through.
11. Add 600 µL of dilution buffer and collect the eluate discarding the first 3 drops (~2 µM).

## A10. Labeling Efficiency

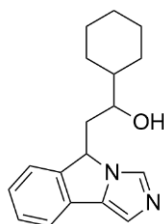
Measurement of protein concentration and degree of labeling (DOL) using a UV-VIS-NIR JASCO V570 Spectrophotometer of JASCO Inc. (Easton, Maryland, USA):

[nanotempertech.com/dol-calculator](https://nanotempertech.com/dol-calculator)

Absorbance A <sub>280</sub>	0.01193	Protein concentration	2.033 µM
Absorbance A <sub>650</sub>	0.01193	Degree-of-labeling (DOL)	1.04

## B1. Ligand/Non-Fluorescent Binding Partner

NLG919 Analogue



## B2. Molecule Class/Organism

Small molecule fragment / IDO1 inhibitor

## B3. Sequence/Formula

$C_{18}H_{22}N_2O$

## B4. Purification Strategy/Source

Selleckchem.com

[S7111](#)

## B5. Stock Concentration/Stock Buffer

14.1 mg/mL | 50 mM

DMSO

## B6. Molecular Weight/Extinction Coefficient

282.4 Da

## B7. Serial Dilution Preparation

1. Add 8  $\mu$ L of DMSO to 2  $\mu$ L of the NLG919 Analogue stock to obtain 10  $\mu$ L of a 10 mM solution.
2. Mix 2  $\mu$ L of the 10 mM NLG919 Analogue solution with 48  $\mu$ L of dilution buffer to obtain 50  $\mu$ L of a 400  $\mu$ M NLG919 Analogue solution.
3. Mix 8  $\mu$ L of DMSO with 192  $\mu$ L of dilution buffer to obtain 200  $\mu$ L of a 4% DMSO solution.
4. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu$ L of the 400  $\mu$ M NLG919 Analogue solution into tube **1**. Then, transfer 10  $\mu$ L of the 4% DMSO solution into tubes **2** to **16**.
5. Prepare a 1:1 serial dilution by transferring 10  $\mu$ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu$ L from tube **16** to get an equal volume of 10  $\mu$ L for all samples.
6. Mix 9  $\mu$ L of labeled IDO1 with 191  $\mu$ L of dilution buffer to obtain 200  $\mu$ L of ~90 nM IDO1.
7. Add 10  $\mu$ L of ~90 nM labeled IDO1 to each tube from **16** to **1** and mix by pipetting.
8. Incubate for 10 minutes at room temperature in the dark before loading capillaries.

## C. Applied Quality Checks

Validation of structural integrity of IDO1 using circular dichroism<sup>1</sup>:

Element	Sample	t = 0'	t = 30'	t = 60'	t = 90'	t = 120'
<b><i>α-helix</i></b>	rhIDO1	28.6%	28.6%	27.2%	27.1%	26.7%
	NT647-rhIDO1	24.5%	24.3%	24.7%	24.0%	24.1%
<b><i>Antiparallel β-sheet</i></b>	rhIDO1	12.7%	12%	13.7%	14.1%	14.5%
	NT647-rhIDO1	15.3%	15.4%	15.0%	15.7%	15.8%
<b><i>Parallel β-sheet</i></b>	rhIDO1	8.9%	8.9%	9.0%	9.0%	9.0%
	NT647-rhIDO1	9.4%	9.4%	9.3%	9.4%	9.3%
<b><i>β-turn</i></b>	rhIDO1	17.5%	17.5%	17.5%	17.5%	17.6%
	NT647-rhIDO1	17.1%	17.2%	17.2%	17.1%	17.2%
<b><i>Random coil</i></b>	rhIDO1	32.3%	32.2%	32.6%	32.3%	32.2%
	NT647-rhIDO1	33.7%	33.7%	33.8%	33.8%	33.6%
<b><i>Total</i></b>	rhIDO1	100%	100%	100%	100%	100%
	NT647-rhIDO1	100%	100%	100%	100%	100%

### D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)

Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH)

### D2. MST Software

MO.Affinity Analysis v2.1 (NanoTemper Technologies GmbH)

### D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

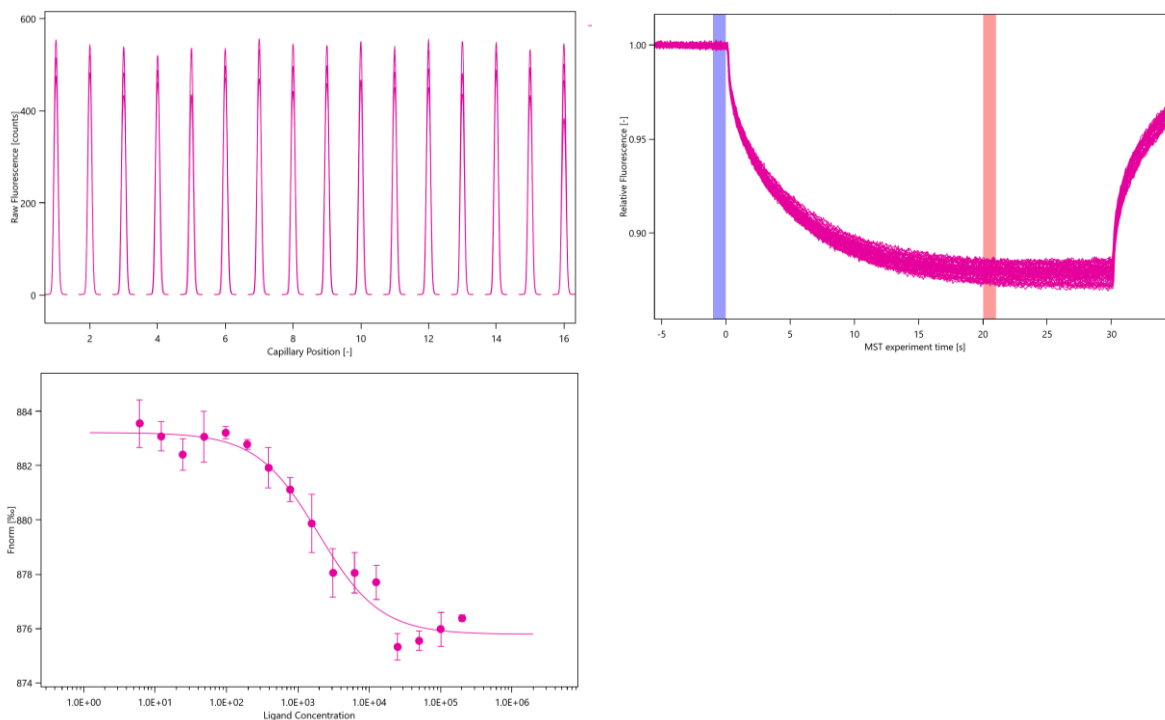
50 mM Tris-HCl, pH 7.4, 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 2 mM DTT, 0.05% TWEEN® 20, 2% DMSO

45 nM IDO1 | 200 μM – 6.1 nM NLG919 Analogue | RT (~23°C) | medium MST power | 20% excitation power

<sup>1</sup> The stability of NT647-rhIDO1 and unmodified rhIDO1 protein was checked using circular dichroism. The spectra of both proteins were recorded using a Jasco810 spectrophotometer with 1 mm path-length quartz cuvettes at room temperature (~22°C). The sensitivity was 100 millidegrees and the scanning speed was 20 nm/min for an accumulation of 2 scans. CD data was collected between 180 and 260 nm for both samples at a concentration of 0.1 mg/mL in phosphate buffer (PPB; 50 mM K<sub>2</sub>HPO<sub>4</sub>, pH 7.4). The deconvolution of the spectra was performed with CDNN 2.1 software.

## D4. MST Results (Capillary Scan/Time Traces/Dose Response)

$K_d = 3.3 \pm 0.4 \mu\text{M}$  [Coletti et al., Eur J Med Chem. 141, 169-177 \(2017\)](#)



## D5. Reference Results/Supporting Results

$EC_{50} = 75 \text{ nM}$  Cell-based activity assay  
[Mautino et al., Proceedings: AACR 104th Annual Meeting 2013; Apr 6-10, 2013](#)

## E. Contributors

Alice Coletti<sup>2</sup>, Antonio Macchiarulo<sup>2</sup>, Semir Jeridi<sup>3</sup>

<sup>2</sup> Università degli Studi di Perugia, Perugia, Italy

<sup>3</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)