

Monolith Protocol MO-P-043

Guanine Quadruplex DNA – 1H6 Monoclonal Antibody

Guanine quadruplex (G4) structures are higher order structures readily formed both *in vitro* and *in vivo* from single stranded guanine-rich RNA or DNA molecules. These structures are often found in the telomeric regions of chromosomes and play a regulatory role in various cellular pathways. 1H6 is a mouse monoclonal antibody which strongly binds to synthetic G4 DNA structures and can serve as a tool for detection of G4 structures *in situ*.

G-quadruplex DNA | antibody | stoichiometry

A1. Target/Fluorescent Molecule

G4 oligonucleotides

A2. Molecule Class/Organism

Synthetic oligonucleotides

A3. Sequence/Formula

(T ₄ G ₄) ₂	5' Cy5 TTT TGG GGT TTT GGG G 3'
(T ₄ G ₃) ₂	5' Cy5 TTT TGG GTT TTG GG 3'
(T ₃ G ₃) ₂	5' Cy5 TTT GGG TTT GGG 3'
(T ₃ G ₄) ₂	5' Cy5 TTT GGG GTT TGG GG 3'
MYC	5' Cy5 TGA GGG TGG GTA GGG TGG GTA A 3'
(A ₄ G ₄) ₂	5' Cy5 AAA AGG GGA AAA GGG G 3'
(C ₄ G ₄) ₂	5' Cy5 CCC CGG GGC CCC GGG G 3'
poly(T)	5' Cy5 TTT TTT TTT TTT TTT 3'

A4. Purification Strategy/Source

Purified with high-performance liquid chromatography (HPLC)
IDT (Leuven, Belgium)

Oligonucleotides were diluted to 10 μM in TE buffer (10 mM Tris-HCl, pH 7.5, 100 mM KCl, 1 mM EDTA). Following denaturation for 10 min at 95°C, G4 structures could form overnight by slowly cooling to room temperature in the heating block.

A5. Stock Concentration/Stock Buffer

100 μM
ddH₂O

A6. Molecular Weight/Extinction Coefficient

5'Cy5	Single stranded, linear (T ₄ G ₄) ₂
532.6 Da	5,005.26 Da
250,000 M ⁻¹ cm ⁻¹ (ε ₆₄₈)	147,000 M ⁻¹ cm ⁻¹ (ε ₂₆₀)

Single stranded sequences are used to prepare quadruplexes. Quadruplexes could be tetramolecular, bimolecular and unimolecular.

A7. Dilution Buffer

400 mM Na PBS, 0.5% BSA, 0.05% TWEEN® 20

A8. Labeling Strategy

5' Cy5 labeled

A9. Labeling Procedure

N/A

A10. Labeling Efficiency

HPLC-purified, 100% labeled DNA

B1. Ligand/Non-Fluorescent Binding Partner

1H6 monoclonal mouse antibody (mAb)

B2. Molecule Class/Organism

Monoclonal antibody

B3. Sequence/Formula

Heavy Chain	CDRH-1				CDRH-2				CDRH-3																	
	27	30	35	38	56	59	62	65	105																	
1H6	G	F	T	F	R	N	Y	W	I	R	L	K	S	D	N	Y	A	T	T	N	W	Y	Y	F	D	Y

Light Chain	CDRL-1				CDRL-2		CDRL-3																
	27	32	34	38	56	65	105																
1H6	Q	S	L	L	Y	S	N	G	K	T	Y	L	V	S	V	Q	G	T	H	F	P	L	T

Adapted from Henderson et al.¹

¹ Henderson et al., *Nucleic Acids Research*, 42 (2), 2014, 860-869

B4. Purification Strategy/Source

Intermolecular Guanine Quadruplex DNA Mouse, Monoclonal Antibody (1H6)
 MédiMabs (Montreal, Canada)
 MM-0265-P

B5. Stock Concentration/Stock Buffer

1 mg/mL | 6.45 μ M
 PBS

B6. Molecular Weight/Extinction Coefficient

150 kDa

B7. Serial Dilution Preparation

1. Mix 31 μ L of the 6.45 μ M 1H6 solution with 69 μ L of dilution buffer to obtain 100 μ L of a 2 μ M solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 40 μ L of the 2 μ M 1H6 solution into tube **1**. Then, transfer 20 μ L of dilution buffer into tubes **2** to **16**.
3. Prepare a 1:1 serial dilution by transferring 20 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 20 μ L from tube **16** to get an equal volume of 20 μ L for all samples.
4. Mix 2 μ L of 10 μ M 5'Cy5 G-quadruplex oligonucleotide with 398 μ L of dilution buffer to obtain 400 μ L of a 50 nM solution.
5. Add 20 μ L of 50 nM 5'Cy5 G-quadruplex oligonucleotide to each tube from **16** to **1** and mix by pipetting.
6. Incubate tubes for 5 minutes at room temperature in the dark before loading capillaries.

Dilution series over wide range to narrow down point of saturation

1. Mix 8 μ L of 10 μ M 5'Cy5 G-quadruplex oligonucleotide with 8 μ L of 10 μ M unlabeled G-quadruplex oligonucleotide and 384 μ L of dilution buffer to obtain 400 μ L of a 400 nM solution.
2. Prepare a PCR-rack with 16 new PCR tubes. Transfer 40 μ L of the 2 μ M 1H6 solution into tube **1**. Then, transfer 20 μ L of dilution buffer into tubes **2** to **16**.
3. Prepare a 1:1 serial dilution by transferring 20 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 20 μ L from tube **16** to get an equal volume of 20 μ L for all samples.
4. Add 20 μ L of 400 nM 5'Cy5 G-quadruplex oligonucleotide to each tube from **16** to **1** and mix by pipetting.
5. Incubate tubes for 5 minutes at room temperature in the dark before loading capillaries.

Dilution series over small concentration range to precisely determine stoichiometry

1. Mix 15 μ L of the 2 μ M 1H6 solution with 123 μ L of dilution buffer to obtain 138 μ L of a 700 nM solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 16 μ L, 15 μ L, ..., 2 μ L, 1 μ L of the 700 nM 1H6 solution into tubes **1** to **16**. Then, transfer 1 μ L, 2 μ L, ..., 14 μ L, 15 μ L of dilution buffer into tubes **2** to **16** to get an equal volume of 16 μ L for all samples.
3. Prepare a 1:1 serial dilution by transferring 16 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 16 μ L from tube **16** to get an equal volume of 16 μ L for all samples.
4. Add 16 μ L of 400 nM 5'Cy5 G-quadruplex oligonucleotide to each tube from **16** to **1** and mix by pipetting.
5. Incubate tubes for 5 minutes at room temperature in the dark before loading capillaries.

D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)
 Capillaries Monolith NT.115 (MO-K022, NanoTemper Technologies GmbH)

D2. MST Software

MO.AffinityAnalysis v2.3 (NanoTemper Technologies GmbH)
nanotempertech.com/monolith/#monolith-software

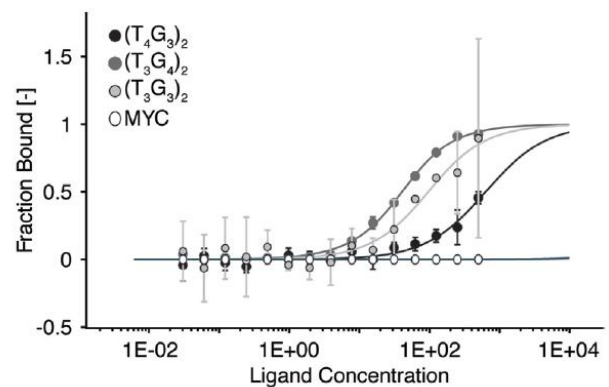
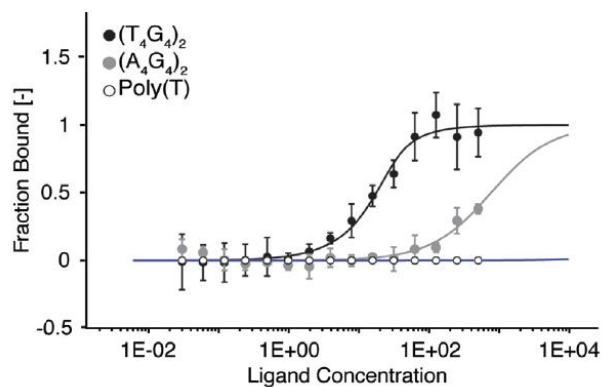
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

400 mM Na PBS, 0.5% BSA, 0.05% TWEEN® 20
 25 nM G4s | 1 µM – 0.03 nM 1H6 antibody | 24°C | medium MST power | 80% excitation power
 400 mM Na PBS, 0.5% BSA, 0.05% TWEEN® 20
 200 nM G4s | 1 µM – 0.03 nM and 350 nM – 50 nM 1H6 antibody | 24°C | medium MST power | 40% excitation power

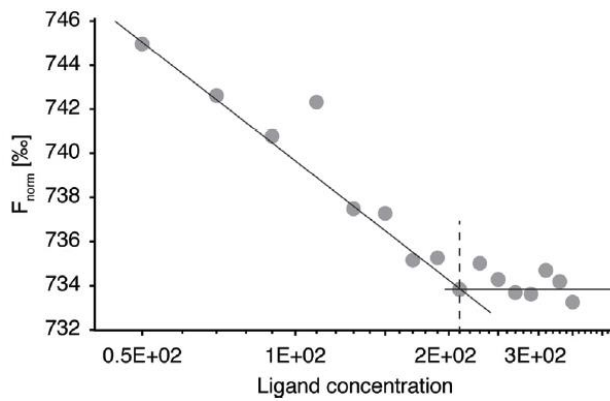
D4. MST Results (Capillary Scan/Time Traces/Dose Response)

$(T_4G_4)_2$	$K_d = 5 \text{ nM}$
$(T_4G_3)_2$	$K_d = 620 \text{ nM}$
$(T_3G_3)_2$	$K_d = 83 \text{ nM}$
$(T_3G_4)_2$	$K_d = 28 \text{ nM}$
MYC	No binding
$(A_4G_4)_2$	$K_d = 754 \text{ nM}$
$(C_4G_4)_2$	$K_d = 184 \text{ nM}$
poly(T)	No binding

Kazemier et al., *Nucleic Acids Research*, 45 (10), 2017, 5913–5919



Intersection of linear fits² at ~215 nM of 1H6 (200 nM of G4S) → 1 : 1 stoichiometry



D5. Reference Results/Supporting Results

G-quadruplex DNA

Sen and Gilbert, *Methods in Enzymology* 211, 1992, 191-199 | Henderson et al., *Nucleic Acids Research*, 42 (2), 2014, 860-869

Hoffmann et al., *Nucleic Acids Research*, 44 (1), 2016, 152-163 | Kazemier et al., *Nucleic Acids Research*, 45 (10), 2017, 5913-5919

E. Contributors

Hinke G. Kazemier³, Katarzyna Walkiewicz⁴

² Linear regression lines of the saturated and non-saturated data points were set manually.

³ European Research Institute for the Biology of Ageing University Medical Center Groningen (ERIBA/UMCG), Netherlands

⁴ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com