

Monolith Protocol MO-P-032

Dendritic Polyglycerols – Human Serum Albumin

Dendritic polyglycerols (dPG) are spherical, water soluble, polyether-based nanomaterials which are decorated with multiple peripheral hydroxyl groups and hold great potential in diagnostics and therapeutic applications. Their high biocompatibility, capability and low toxicity *in vitro* as well as *in vivo* make them attractive nanomaterials in the field of nanotheranostics. To evaluate their potential, it is important to understand their cellular behavior and physiological fate (e.g. absorption, metabolism, distribution and excretion). The interaction with serum proteins such as human serum albumin is therefore an important part in the analysis of dPG.

protein – small molecule interaction | theranostics

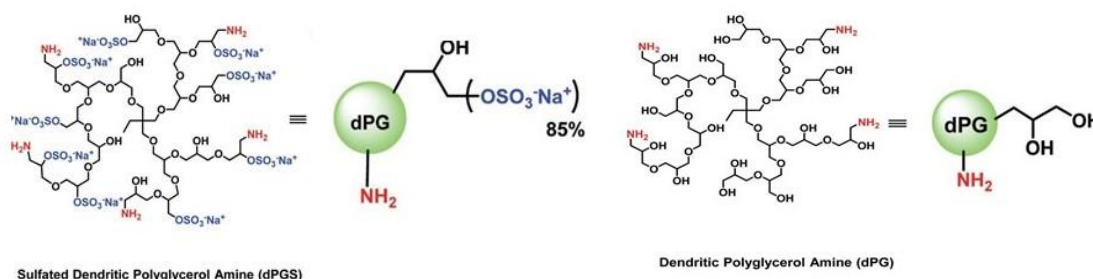
A1. Target/Fluorescent Molecule

Sulfated dPG (dPGS)

A2. Molecule Class/Organism

Dendritic polyglycerol (dPG)

A3. Sequence/Formula



A4. Purification Strategy/Source

Preparation of dendritic polyglycerol (dPG) and their corresponding sulfated (dPGS) amines according to procedures previously reported.¹

A5. Stock Concentration/Stock Buffer

N/A

A6. Molecular Weight/Extinction Coefficient

22 kDa (dPGS) / 10 kDa (dPG)

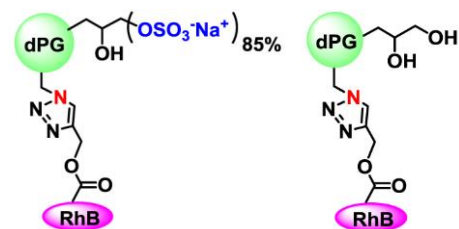
¹ Sunder et al., *Macromolecules* 32 (1999) 4240–4246 | Frey et al., *Mol Biotechnol* 90 (2002) 257–267 | Weinhart et al., *Macromolecules* 11 (2011) 1088–1098

A7. Dilution Buffer

Phosphate buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

A8. Labeling Strategy

Conjugation of a rhodamine B (RhB) dye to azide terminated dPGS/dPG via click chemistry



dPGS-RhB (5b)

dPG(OH)-RhB (2b), negative control

A9. Labeling Procedure

RhB was conjugated to dPGS/dPG according to a previously published procedure.²

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Human serum albumin (HSA)

uniprot.org/uniprot/P02768

B2. Molecule Class/Organism

Serum protein

Homo sapiens (Human)

B3. Sequence/Formula

DAHKSEVAHR FKDLGEENFK ALVLIAFAQY LQQCPFEDHV KLVNEVTEFA KTCVADESAE NCDKSLHTLF GDKLCTVATL
 RETYGEMADC CAKQEPERNE CFLQHKDDNP NLPRLVRPEV DVMCTAFHDN EETFLKQYLY EIARRHPYFY APELLFFAKR
 YKAAFTTECCQ AADKAACLLP KLDEL RDEGK ASSAKQRLKC ASLQKFGERA FKAWAVARLS QRFPAEFAE VSKLVDLTK
 VHTECCHGDL LECADDRADL AKYICENQDS ISSKLKECE KPLLEKSHCI AEVENDEMPA DLPSLAADFV ESKDVCKNYA
 EAKDVF LGMF LYEYARRHPD YSVVLLRLA KTYETTLEKC CAAADPHECY AKVFDEFKPL VEEPQNLIKQ NCELFEQLGE
 YKFQNALVR YTKKVPQVST PTLVEVSRNL GKVGSKCCKH PEAKRMPCAE DYLSVVLNQL CVLHEKTPVS DRVTKCCTES
 LVNRRPCFSA LEVDETYVPK EFNAETTFH ADICTLSEKE RQIKKQTALV ELVKHKPKAT KEQLKAVMDD FAAFVEKCKC
 ADDKETCF AE EGKLV AASQ AALGL

² Gröger et al., *Adv Healthcare Mater* 3 (2014) 375-385

B4. Purification Strategy/Source

Sigma-Aldrich GmbH

SRP6182

B5. Stock Concentration/Stock Buffer

500 μM

Phosphate buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

B6. Molecular Weight/Extinction Coefficient

66.5 kDa

43,800 $\text{M}^{-1}\text{cm}^{-1}$ (ϵ_{280})

B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 500 μM HSA solution into tube **1**. Then, transfer 10 μL of dilution buffer into tubes **2** to **16**.
 2. Prepare a 1:1 serial dilution by transferring 10 μL from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μL from tube **16** to get an equal volume of 10 μL for all samples.
 3. Add 10 μL of 30 nM dPG(S)-RhB or dPG(OH)-RhB to each tube from **16** to **1** and mix by pipetting.
 4. Incubate 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.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

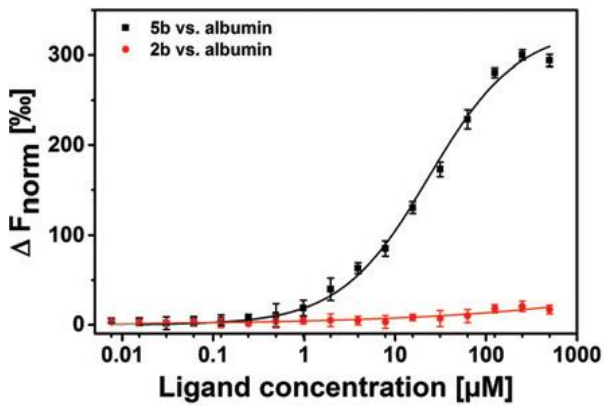
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

15 nM dPGS | 250 μM – 7.6 nM HSA | 25°C

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

EC₅₀ = 12 μM Pant et al., *Nanoscale* 9 (2017) 8723–8739



Interaction of HSA with dPG(S)-RhB (5b, black) and dPG(OH)-RhB (2b, red, negative control).

D5. Reference Results/Supporting Results

N/A

E. Contributors

Kritee Pant³, Teresia Hallström⁴

³ Institute of Radiopharmaceutical Cancer Research, Helmholtz-Zentrum, Dresden, Germany

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