



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

CALXGLUK

(Cat. # 786-1524, 786-1525 & 786-1526)



think proteins! think G-Biosciences www.GBiosciences.com

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1524	CALXGLUK	10 mg
786-1525	CALXGLUK	50 mg
786-1526	CALXGLUK	100mg

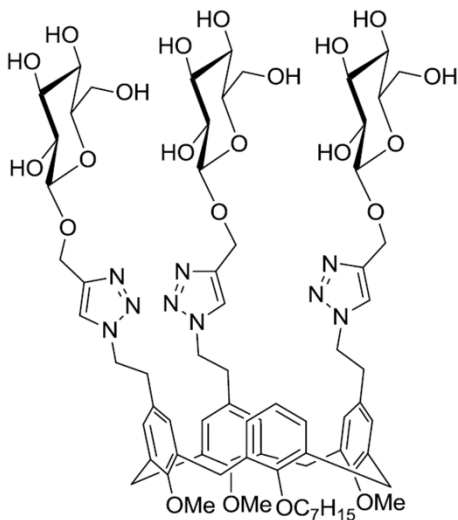
REAGENT INFORMATION

- Compound Name: CALXGLUK
- IUPAC Name: 5,11,17-tris{2-[4-(β-D-glucopyranosyl)-1H-1,2,3-triazol-1-yl]ethyl}-25-monoheptoxy-26,27,28-trimethoxycalix[4]arene
- Application: Membrane Protein Extraction
- Surfactant type: Calixarene
- Molecular Formula: C₇₁H₉₅N₉O₂₂
- Molecular Weight: 1426.5 g/mol
- Percent composition: C (59.78%), H (6.71%), N (8.84%), O (24.67%)
- Physical state: Beige powder
- Purity (HPLC, 280nm): <95%
- Retention Time (RP₁₈ HPLC): t_R = 13.4 min
- Critical Micelle Concentration: 0.025mM

STORAGE AND HANDLING

- Storage conditions: Store in <-20°C freezer for up to one year
- Solubility: Soluble in water (2.5mM), methanol and DMSO.

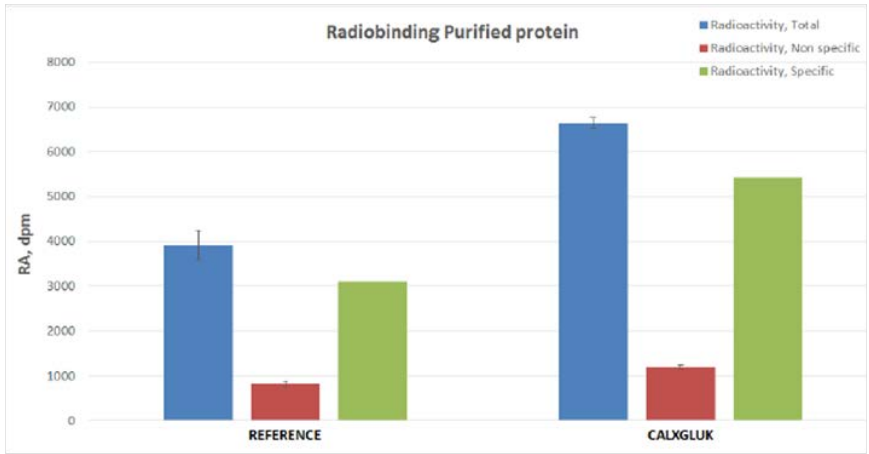
STRUCTURE



BIOCHEMICAL VALIDATION DATA

Binding of radioligand on target protein, purified in reference detergent with or without CALXGLUK as an additive.

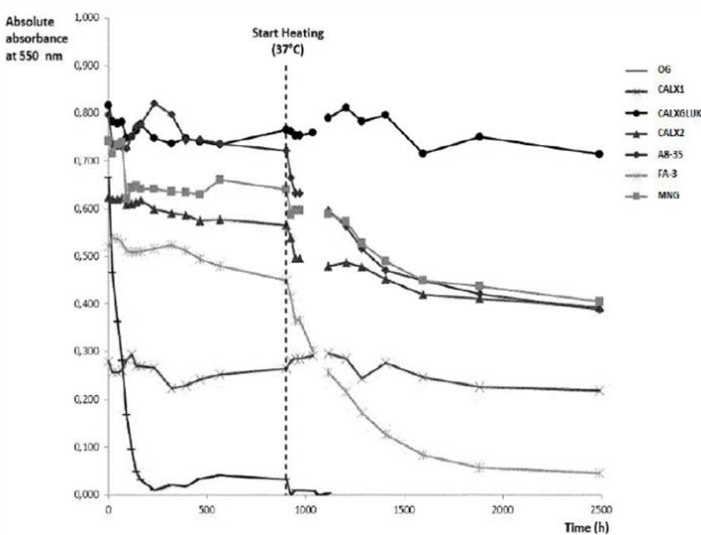
Purified protein was incubated with radioligand in absence (total, blue bars) or presence (Non Specific signal, red bars) of an excess of cold ligand. After filtration on GF/C membranes and washing, scintillation agent was added and radioactivity was detected using a Microbeta2. Specific radioactivity (green bars) corresponds to (total signal) – (non-specific signal).



Stabilization of the Bacteriorhodopsin

After solubilization and purification in OG, reagents were added as stabilizers.

Bacteriorhodopsin's activity was monitored over time at first 25°C and then 37°C by measuring absorbance at 560nm.



REFERENCES

1. Proprietary patent: "Method for the synthesis of steroidal calixarene compounds and their use for extraction, solubilization and/or stabilization of native and functional membrane proteins"; WO 2017134087 A1 ; Publication date: August 10, 2017.

Last saved: 7/17/2018 EM



www.GBiosciences.com