



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

A Geno Technology, Inc. (USA) brand name

FTAC6

(Cat. # 786-1530, 786-1531 & 786-1532)



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

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1530	FTAC6	250 mg
786-1531	FTAC6	500 mg
786-1532	FTAC6	1 g

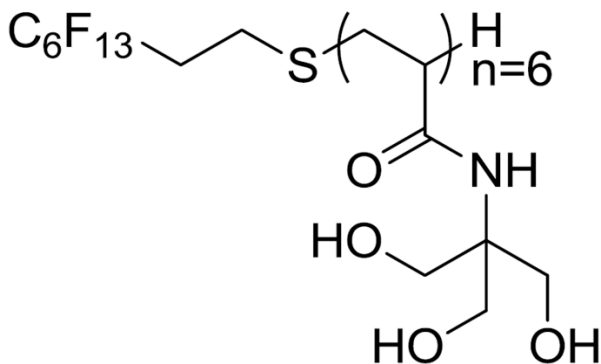
REAGENT INFORMATION

- Compound Name: FTAC6
- IUPAC Name: S-(poly(tris(hydroxymethyl)acrylamidomethane)-(3,3,4,4,5,5,6,6,7,7,8,8,8-tridecafluorooctyl) DP_n=6
- Application: Membrane Protein Extraction
- Molecular Formula: na
- Molecular Weight: ≈1400 g/mol
- Physical state: White powder
- Retention Time (RP₁₈ HPLC): t_R = 10.5 min
- Critical Micelle Concentration: 0.37mM

STORAGE AND HANDLING

- Storage conditions: Store in <-20°C freezer for up to one year
- Solubility: Soluble in water (18.5mM), methanol and DMSO.

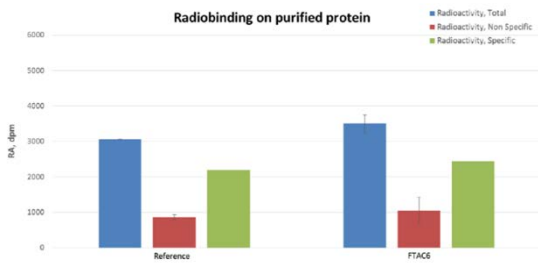
STRUCTURE



BIOCHEMICAL VALIDATION DATA

Binding of radioligand on GPCR protein, purified in reference detergent with or without addition of FTAC6 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)



REFERENCES

1. Damian, M., S. Perino, A. Polidori, A. Martin, L. Serre, B. Pucci and J.-L. Baneres (2007). "New tensio-active molecules stabilize a human G protein-coupled receptor in solution." *Febs Letters* 581(10): 1944-1950.
2. Park, K.-H., C. Berrier, F. Lebaupain, B. Pucci, J.-L. Popot, A. Ghazi and F. Zito (2007). "Fluorinated and hemifluorinated surfactants as alternatives to detergents for membrane protein cell-free synthesis." *Biochemical Journal* 403: 183-187.
3. Joubert, O., R. Nehme, D. Fleury, M. De Rivoyre, M. Bidet, A. Polidori, M. Ruat, B. Pucci, P. Mollat and I. Mus-Veteau (2009). "Functional studies of membranebound and purified human Hedgehog receptor Patched expressed in yeast." *Biochimica Et Biophysica Acta-Biomembranes* 1788(9): 1813-1821.
4. Nehme, R., O. Joubert, M. Bidet, B. Lacombe, A. Polidori, B. Pucci and I. MusVeteau (2010). "Stability study of the human G-protein coupled receptor, Smoothened." *Biochimica Et Biophysica Acta-Biomembranes* 1798(6): 1100-1110.
5. Park, K.-H., E. Billon-Denis, T. Dahmane, F. Lebaupain, B. Pucci, C. Breyton and F. Zito (2011). "In the cauldron of cell-free synthesis of membrane proteins: playing with new surfactants." *New Biotechnology* 28(3): 255-261.
6. Kyrychenko, A., M. V. Rodnin, M. Vargas-Urbe, S. K. Sharma, G. Durand, B. Pucci, J.-L. Popot and A. S. Ladokhin (2012). "Folding of diphtheria toxin T-domain in the presence of amphipols and fluorinated surfactants: Toward thermodynamic measurements of membrane protein folding." *Biochimica Et Biophysica ActaBiomembranes* 1818(4): 1006-1012.

Last saved: 7/17/2018 EM



www.GBiosciences.com