

PRODUCT INFORMATION



FTAC6

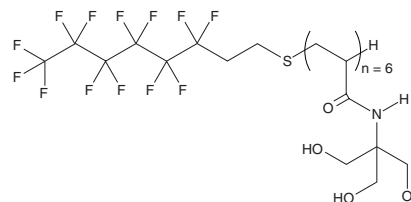
Item No. 24786

Formal Name: N¹,N³,N⁵,N⁷,N⁹,N¹¹-hexakis(1,3-dihydroxy-2-(hydroxymethyl)propan-2-yl)-12-((3,3,4,4,5,5,6,6,7,7,8,8,8-tridecafluorooctyl)thio)dodecane-1,3,5,7,9,11-hexacarboxamide

Supplied as: A powder

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

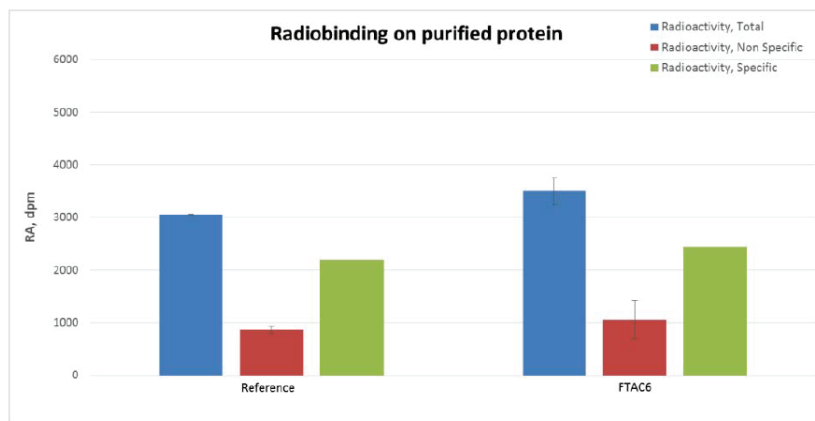
Laboratory Procedures

FTAC6 is supplied as a powder. A stock solution may be made by dissolving the FTAC6 in the solvent of choice, which should be purged with an inert gas. FTAC6 is soluble in organic solvents such as methanol and DMSO. FTAC6 is also soluble in water at a concentration of approximately 18.5 mM. We do not recommend storing the aqueous solution for more than one day.

Description

FTAC6 is a detergent that can be used to stabilize membrane proteins. It has a critical micelle concentration (CMC) of 0.37 mM. It has been used in the cell-free synthesis and purification of histidine-tagged MscL, a mechanosensitive membrane channel, as well as in the insertion of MscL into liposomes.¹ FTAC6 has also been used in the synthesis of MscL *in vitro* and preserves its activity.

Image



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).

Reference

1. Park, K.-H., Berrier, C., Lebaupain, F., *et al.* Fluorinated and hemifluorinated surfactants as alternatives to detergents for membrane protein cell-free synthesis. *Biochem. J.* **403**(1), 183-187 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM