

# PRODUCT INFORMATION

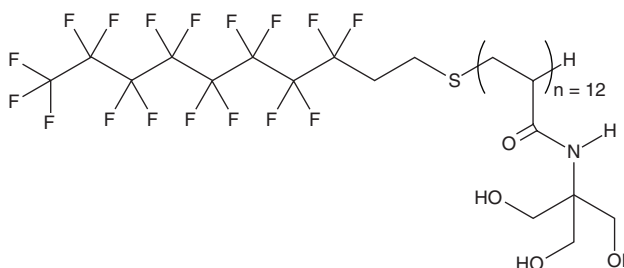


## FTAC8

Item No. 24787

**Formal Name:** N<sup>1</sup>,N<sup>3</sup>,N<sup>5</sup>,N<sup>7</sup>,N<sup>9</sup>,N<sup>11</sup>,N<sup>13</sup>,N<sup>15</sup>,N<sup>17</sup>,N<sup>19</sup>,N<sup>21</sup>,N<sup>23</sup>-dodecakis(1,3-dihydroxy-2-(hydroxymethyl)propan-2-yl)-24-((3,3,4,4,5,5,6,6,7,7,8,8,9,9,10,10,10-heptadecafluorodecyl)thio)tetracosane-1,3,5,7,9,11,13,15,17,19,21,23-dodecacarboxamide

**Synonym:** C<sub>8</sub>F-TAC  
**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.

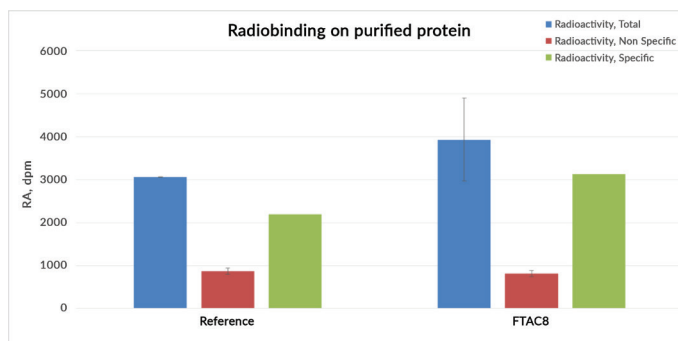
## Laboratory Procedures

FTAC8 is supplied as a powder. A stock solution may be made by dissolving the FTAC8 in the solvent of choice, which should be purged with an inert gas. FTAC8 is soluble in organic solvents such as methanol and DMSO. It is also soluble in water at a concentration of 1 mM. We do not recommend storing the aqueous solution for more than one day.

## Description

FTAC8 is a detergent that can be used to stabilize membrane proteins. It has a critical micelle concentration (CMC) of 0.02 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.<sup>1</sup> FTAC8 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 FTAC8 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