

PRODUCT INFORMATION



FLAC6

Item No. 24781

CAS Registry No.: 2820046-26-0
Formal Name: (2R,3R,4R,5R)-2,3,5,6-tetrahydroxy-N-(4,4,5,5,6,6,7,7,8,8,9,9,9-tridecafluorononyl)-4-(((2S,3R,4S,5R,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)hexanamide

MF: C₂₁H₂₈F₁₃NO₁₁

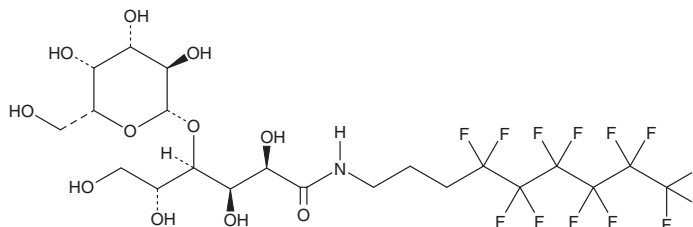
FW: 717.4

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

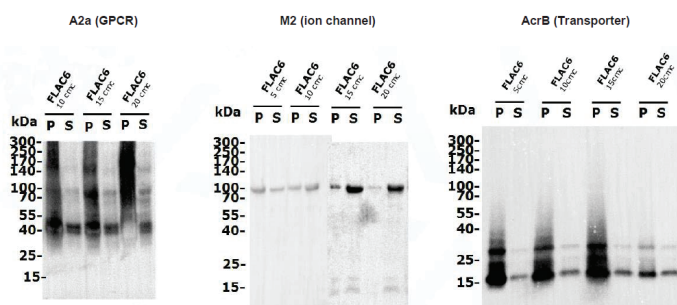
FLAC6 is supplied as a crystalline solid. A stock solution may be made by dissolving the FLAC6 in the solvent of choice, which should be purged with an inert gas. FLAC6 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of FLAC6 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of FLAC6 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

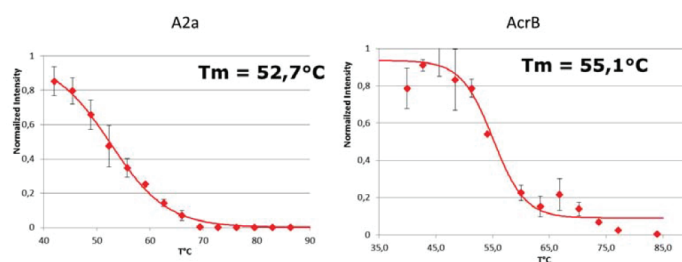
Description

FLAC6 is a detergent that can be used to solubilize membrane proteins. It has a critical micelle concentration (CMC) of 0.56 mM.

Images



Membrane protein solubilization from various membranes. Various targets were extracted from biological membranes by using FLAC6 reagent at 5- to 20-fold the critical micelle concentration (CMC). After solubilization, samples were centrifuged at 100,000 g. Proteins from pellets (P) and supernatants (S) were separated on a 4-15% Tris-glycine SDS-PAGE, transferred to PVDF membrane and immunodetected with either specific or anti-tag antibodies.
T=total, P=pellet, S=supernatant



Thermal assay on membranes solubilized with 15-fold FLAC6 CMC. Unfolding of either A2a receptor or AcrB was followed by Western blotting after applying a temperature gradient and high speed centrifugation.

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, 01/06/2023

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