

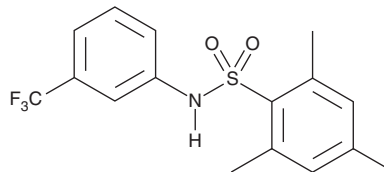
PRODUCT INFORMATION



m-3M3FBS

Item No. 16867

CAS Registry No.: 200933-14-8
Formal Name: 2,4,6-trimethyl-N-[3-(trifluoromethyl)phenyl]-benzenesulfonamide
MF: C₁₆H₁₆F₃NO₂S
FW: 343.4
Purity: ≥95%
UV/Vis.: λ_{max}: 230 nm
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

m-3M3FBS is supplied as a crystalline solid. A stock solution may be made by dissolving the *m*-3M3FBS in the solvent of choice, which should be purged with an inert gas. *m*-3M3FBS is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *m*-3M3FBS in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

m-3M3FBS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, *m*-3M3FBS should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. *m*-3M3FBS has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

m-3M3FBS is an activator of phospholipase C (PLC) that stimulates superoxide generation, increase in cytoplasmic calcium, and inositol phosphate formation in human neutrophils when used at a concentration of 15-50 μM.¹ It stimulates cytoplasmic calcium increases in several cell lines and, at 25 μM *in vitro*, promotes the hydrolysis of phosphatidylinositol bisphosphate by all PLC isoforms.¹ *m*-3M3FBS is used to study PLC signaling in cells and animals, often in conjunction with the PLC inhibitor U-73122 (Item No. 70740).²⁻⁴

References

1. Bae, Y.S., Lee, T.G., Park, J.C., *et al.* Identification of a compound that directly stimulates phospholipase C activity. *Mol. Pharmacol.* **63**(5), 1043-1050 (2003).
2. Kim, S.D., Kim, H.J., Shim, J.W., *et al.* Phospholipase C activator *m*-3M3FBS protects against morbidity and mortality associated with sepsis. *J. Immunol.* **189**(4), 2000-2005 (2012).
3. Semple-Rowland, S., Madorsky, I., Bolch, S., *et al.* Activation of phospholipase C mimics the phase shifting effects of light on melatonin rhythms in retinal photoreceptors. *PLoS One* **8**(12), e83378 (2013).
4. Szebenyi, S.A., Ogura, T., Sathyanesan, A., *et al.* Increases in intracellular calcium via activation of potentially multiple phospholipase C isozymes in mouse olfactory neurons. *Front. Cell. Neurosci.* **8**, 336 (2014).

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, 09/30/2022

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