

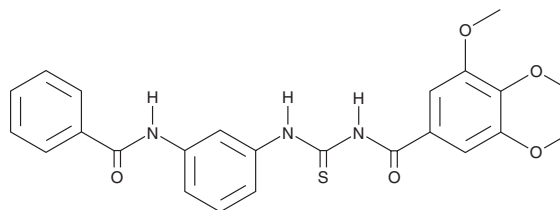
PRODUCT INFORMATION



MRT10

Item No. 22138

CAS Registry No.: 330829-30-6
Formal Name: N-[[[3-(benzoylamino)phenyl]amino]thioxomethyl]-3,4,5-trimethoxy-benzamide
MF: C₂₄H₂₃N₃O₅S
FW: 465.5
Purity: ≥98%
UV/Vis.: λ_{max}: 284 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

MRT10 is supplied as a crystalline solid. A stock solution may be made by dissolving the MRT10 in the solvent of choice, which should be purged with an inert gas. MRT10 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MRT10 in these solvents is approximately 33 mg/ml.

MRT10 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MRT10 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MRT10 has a solubility of approximately 0.3 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MRT10 is an antagonist of the Smoothed (Smo) receptor with an IC₅₀ value of 0.64 μM in a luciferase reporter assay.^{1,2} It also inhibits bopipy-cyclopamine binding to the murine Smo receptor (IC₅₀ = 0.5 μM) when expressed in HEK293 cells.¹

References

1. Manetti, F., Faure, H., Roudaut, H., *et al.* Virtual screening-based discovery and mechanistic characterization of the acylthiourea MRT-10 family as smoothed antagonists. *Mol. Pharm.* **78(4)**, 658-665 (2010).
2. Solinas, A., Faure, H., Roudaut, H., *et al.* Acylthiourea, acylurea, and acylguanidine derivatives with potent hedgehog inhibiting activity. *J. Med. Chem.* **55(4)**, 1559-1571 (2012).

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, 11/07/2022

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