

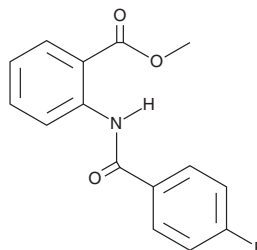
PRODUCT INFORMATION



Exo1

Item No. 18455

CAS Registry No.: 75541-83-2
Formal Name: 2-[(4-fluorobenzoyl)amino]-benzoic acid, methyl ester
Synonyms: Exocytic Inhibitor 1, NSC 214045
MF: C₁₅H₁₂FNO₃
FW: 273.3
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 267, 314 nm
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Exo1 is supplied as a solid. A stock solution may be made by dissolving the Exo1 in the solvent of choice, which should be purged with an inert gas. Exo1 is soluble in organic solvents such as ethanol and DMSO. The solubility of Exo1 in these solvents is approximately 25 and 100 mM, respectively.

Exo1 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Exo1 is a cell-permeable, reversible inhibitor of exocytosis (IC₅₀ = 20 μM).¹ It interferes with protein trafficking and secretion mediated by the Golgi apparatus and endoplasmic reticulum by dissociating the GTPase activator ADP-ribosylation factor (ARF1) from Golgi membranes, which leads to a consequent collapse of the Golgi apparatus.¹

Reference

1. Feng, Y., Yu, S., Lasell, T.K.R., et al. Exo1: A new chemical inhibitor of the exocytic pathway. *Proc. Natl. Acad. Sci. USA* **100(11)**, 6469-6474 (2003).

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, 10/11/2022

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