

PRODUCT INFORMATION

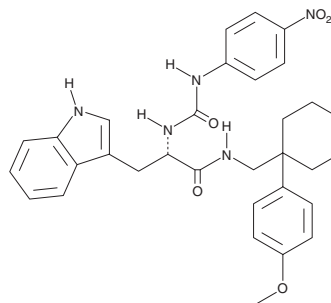


ML-18

Item No. 27319

CAS Registry No.: 1422269-30-4
Formal Name: (αS)-N-[[1-(4-methoxyphenyl)cyclohexyl)methyl]-α-[[[(4-nitrophenyl)amino]carbonyl]amino]-1H-indole-3-propanamide

MF: C₃₂H₃₅N₅O₅
FW: 569.7
Purity: ≥98%
Supplied as: A 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

ML-18 is supplied as a solid. A stock solution may be made by dissolving the ML-18 in the solvent of choice, which should be purged with an inert gas. ML-18 is soluble in the organic solvent DMSO.

Description

ML-18 is a non-peptide bombesin receptor subtype 3 (BB₃) antagonist that inhibits [¹²⁵I]BA1 binding to NCI-H1299 lung cancer cells transfected with human BB₃ receptors (IC₅₀ = 4.8 μM).¹ It is selective for BB₃ over the gastrin releasing peptide receptor (GRPR) and the neuromedin B receptor (NMBR) with IC₅₀ values of 16 and >100 μM, respectively, in a radioligand binding assay. ML-18 (16 μM) reversibly inhibits BA1-induced increases in cytosolic calcium in NCI-H1299 cells. It also inhibits BA1-induced phosphorylation of ERK and EGFR and reduces proliferation of NCI-H1299 cells.

Reference

1. Moody, T.W., Mantey, S.A., Moreno, P., *et al.* ML-18 is a non-peptide bombesin receptor subtype-3 antagonist which inhibits lung cancer growth. *Peptides* 64, 55-61 (2015).

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, 11/30/2022

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