

Exo1

ADP-ribosylation factor inhibitor

Exo1 inhibits traffic emanating from the endoplasmic reticulum by inducing rapid collapse of the Golgi to the endoplasmic reticulum¹. Like Brefeldin A, Exo1 induces the rapid release of ARF1 from Golgi membranes, but evidence suggests that Exo1 acts by a different mechanism. Exo1 does not induce the ADP-ribosylation of CtBP/Bars50, and does not interfere with the activity of guanine nucleotide exchange factors specific for Golgi-based ARFs, allowing Exo1 to be used as a tool for differentiating the fatty acid exchange activity of Bars50 from ARF1 activity.

Citations: 1

[View Online »](#)

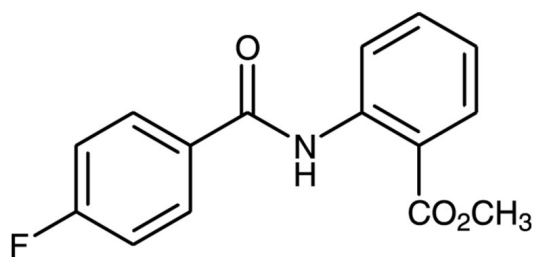
Ordering Information

[Order Online »](#)

BML-EI322-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-[(4-Fluorobenzoyl)amino]benzoic acid methyl ester
Appearance	White solid.
CAS	461681-88-9
Couple Target	ADP-ribosylation factor
Couple Type	Inhibitor
Formula	$C_{15}H_{12}FNO_3$
MW	273.3
MeltingPoint	132-134°C
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10 mg/ml warm), DMF (20 mg/ml), or ethanol (5 mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)