

CCG-1423

Rho inhibitor

Cell permeable inhibitor of RhoA- and RhoC-mediated signaling. Inhibits LPA-induced DNA synthesis in PC-3 prostate cancer cell lines (<1 μ M) and inhibits the growth of RhoC-overexpressing melanoma cell lines (A375M2 and SK-Mel-147) at nanomolar concentrations.

Citations: 2

[View Online »](#)

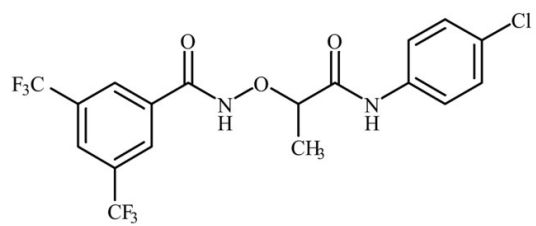
Ordering Information

[Order Online »](#)

BML-EI394-0010	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	N-(2-(4-Chloroanilino)-1-methyl-2-oxoethoxy)-3,5-bis(trifluoromethyl)benzamide
Appearance	White solid.
CAS	285986-88-1
Couple Target	Rho
Couple Type	Inhibitor
Formula	$C_{18}H_{13}ClF_6N_2O_3$
MW	454.8
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol (9mg/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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)