

DIF-1

Ultra-pure; Induces cell differentiation

Supresses cell growth and promotes retinoic acid-induced cell differentiation in HL-60 cells. It is a known inhibitor of Wnt/ β -Catenin signaling pathway.

Citations: 4

[View Online »](#)

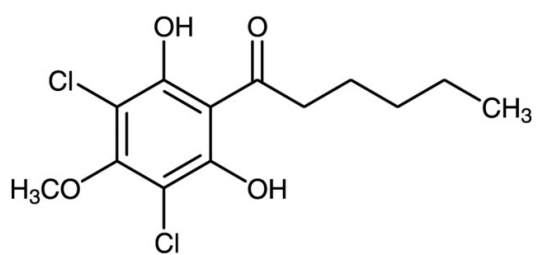
Ordering Information

[Order Online »](#)

BML-GR324-0100	100 μ g
BML-GR324-0500	500 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-(3,5-Dichloro-2,6-dihydroxy-4-methoxyphenyl) hexan-1-one, Differentiation-inducing factor-1

Appearance Pale yellow solid.

CAS 111050-72-7

Couple Target Wnt

Couple Type Inhibitor

Formula $C_{13}H_{16}Cl_2O_4$

MW 307.2

Purity ≥98% (HPLC)

Solubility Soluble in ethanol.

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)