

L-858051

Adenylate cyclase activator

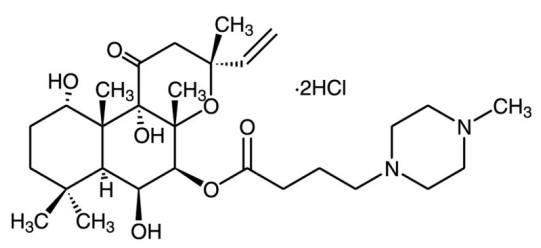
A water-soluble forskolin derivative, retaining adenylate cyclase stimulatory activity ($EC_{50}=10 \mu\text{M}$).

Citations: 5

[View Online »](#)

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

Couple Target Adenylate cyclase

Couple Type Activator

Formula $C_{29}H_{48}N_2O_7 \cdot 2HCl \cdot H_2O$

MW 627.6

Purity ≥92% (HPLC)

Solubility Soluble in water (>10 mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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