

Diacylglycerol Kinase Inhibitor II (R59 949)

Potentiates PKC activity

Inhibitor of diacylglycerol kinase in intact cells ($IC_{50}=0.12\mu M$) and in isolated membranes ($IC_{50}=0.3\mu M$). More potent analog of R59 022 (Prod. No. ALX-430-027). Useful for elucidating roles of protein kinase C (PKC).

Citations: 5

[View Online »](#)

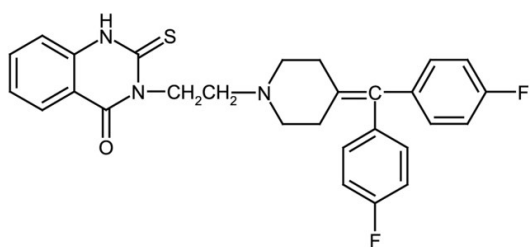
Ordering Information

[Order Online »](#)

ALX-430-028-M005	5mg
ALX-430-028-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3-[2-[4-[bis(4-Fluorophenyl)methylene]-1-piperidinyl]ethyl]-2,3-dihydro-2-thioxo-4(1H)-quinazolinone

Appearance White to off-white powder.

CAS 120166-69-0

Couple Target Diacylglycerol kinase

Couple Type Inhibitor

Formula $C_{28}H_{25}F_2N_3OS$

MW 489.6

Purity ≥98% (HPLC)

Solubility Soluble in dichloromethane, ethyl acetate, tetrahydrofuran, DMSO (33mg/ml) or dimethyl formamide; almost insoluble in water.

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-be@enzolifesciences.com

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