

Gö 6976

PKC inhibitor

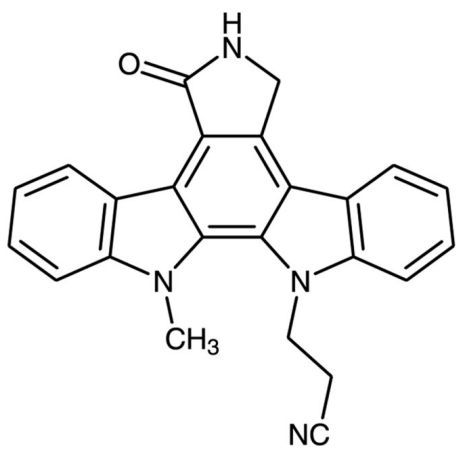
Very potent protein kinase C (PKC) inhibitor with high selectivity for the PKC isotypes α and β . Inhibits PKC α (IC_{50} =2.3nM) and PKC β 1 (IC_{50} =6.2nM) but has no effect on δ , ϵ or γ isotypes. Potent antagonist of HIV-1 induction.

Citations: 8

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for at least 3-4 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	12-(2-Cyanoethyl)-6,7,12,13-tetrahydro-13-methyl-5-oxo-5H-indolo[2,3-a]pyrrolo[3,4-c]carbazole
Appearance	Off-white to yellow solid.
CAS	136194-77-9
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{24}H_{18}N_4O$
MW	378.4
Purity	≥95% (HPLC, TLC)
Solubility	Soluble in DMSO (25mg/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