

PKC-412

Flt inhibitor

Inhibitor of a variety of serine/threonine and tyrosine kinases, like protein kinase C (PKC), cyclic AMP-dependent protein kinase (PKA), S6 kinase, Akt (protein kinase B; PKB), epidermal growth factor receptor (EGFR) tyrosine kinase activity and others including KDR, VEGFR, PDGFR, c-kit and other receptor tyrosine kinases. Potently inhibits FLT-3 kinase including mutant forms found in acute myeloid leukemia *in vitro* and *in vivo*. Apoptosis inducer. Showed broad antiproliferative activity against various tumor cell lines. Selectively inhibits T lymphocyte production of TNF- α . Upregulates endothelial nitric oxide synthase (eNOS; NOS III). Abrogates tumor angiogenesis *in vivo*.

Citations: 29

[View Online »](#)

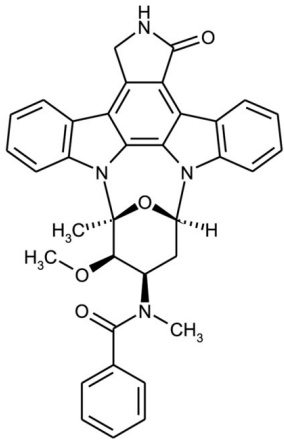
Ordering Information

[Order Online »](#)

BML-EI370-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4'-N-Benzoyl-staurosporine, Midostaurin, CGP 41251
Appearance	White to off-white powder.
CAS	120685-11-2
Couple Target	Akt, c-Kit, EGFR, Flt, PDGF receptor, PKA, PKC, S6 kinase, VEGFR
Couple Type	Inhibitor
Formula	$C_{35}H_{30}N_4O_4$
MI	14: 6185
MW	570.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (15mg/ml), MDC (10mg/ml), methanol (5mg/ml) or 100% ethanol (2.5mg/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-
be@enzolifesciences.com](mailto:info-be@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)