

Erbstatin Analog

EGFR inhibitor

Erbstatin is a novel inhibitor of the EGF receptor-associated tyrosine kinase. It is an unstable compound and is completely inactivated in serum in 30 minutes. Methyl 2,5-dihydroxycinnamate is a stable erbstatin analog retaining activity after a 60 minute incubation. It inhibits EGF receptor-associated tyrosine kinase *in vitro* ($IC_{50}=0.77\text{ }\mu\text{M}$). Inhibition is competitive with substrate and noncompetitive with ATP.

Citations: 5

[View Online »](#)

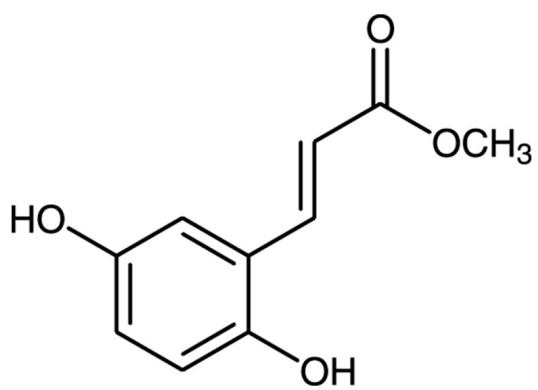
Ordering Information

[Order Online »](#)

BML-EI146-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light and oxygen.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Methyl 2,5-dihydroxycinnamate
Appearance	Yellow solid.
CAS	63177-57-1
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{10}H_{10}O_4$
MW	194.2
MeltingPoint	169-172°C
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (25mg/ml), methanol or DMSO (25mg/ml).



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)