

# 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 »](#)

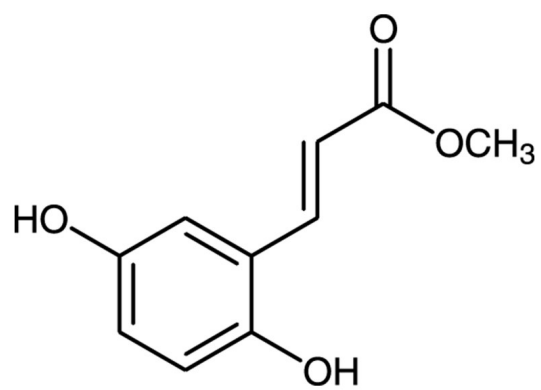
## Ordering Information

[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). |

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)