

Tyrphostin 23 (RG-50810)

EGFR inhibitor

A potent inhibitor of EGF receptor kinase activity ($IC_{50}=35\text{ }\mu\text{M}$).

Citations: 6

[View Online »](#)

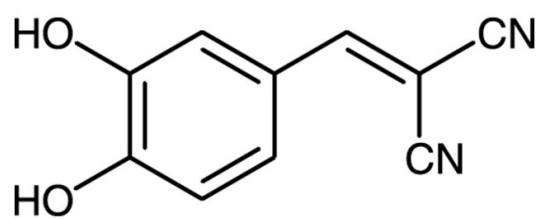
Ordering Information

[Order Online »](#)

BML-EI191-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Product is subject to oxidation.
Handling	Store tightly sealed in the dark. Protect from light. Make solutions fresh daily.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	α -Cyano-(3,4-dihydroxy)cinnamonnitrile
Appearance	Yellow-tan solid.
CAS	118409-57-7
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{10}H_6N_2O_2$
MW	186.1
Purity	≥99% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)