

Tyrphostin 8

Calcineurin inhibitor

The tyrphostins are well known tyrosine kinase inhibitors. However, tyrphostin 8 has been shown to inhibit calcineurin ($IC_{50}=21\text{ }\mu\text{M}$) while displaying weak inhibition of EGFR tyrosine kinase ($IC_{50}=560\text{ }\mu\text{M}$).

Citations: 4

[View Online »](#)

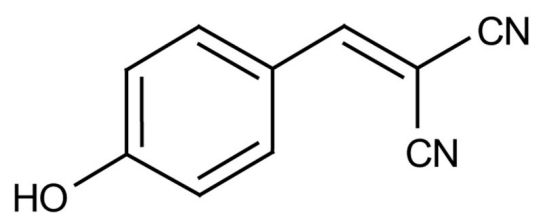
Ordering Information

[Order Online »](#)

BML-PR113-0010	10mg
BML-PR113-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Store tightly sealed in the dark. Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	α -Cyano-(4-hydroxy)cinnamitrile
Appearance	Light orange solid.
CAS	3785-90-8
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{10}H_6N_2O$
MW	170.2
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (25mg/ml) 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)