

Tyrphostin AG 527

Citations: 1

[View Online »](#)

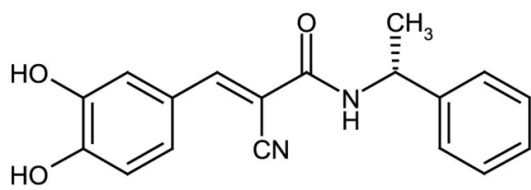
Ordering Information

[Order Online »](#)

ALX-270-141-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Store tightly sealed in the dark. Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tyrphostin B44, (-)-(R)-N-(α -Methylbenzyl)-3,4-dihydroxybenzylidenecyanoacetamide

Appearance Yellow solid.

CAS 133550-32-0

Formula $C_{18}H_{16}N_2O_3$

MW 308.3

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in DMSO or 100% ethanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com