

Tomatidine . HCl

Negative control for Cyclopamine

Steroidal alkaloid structurally similar to cyclopamine (Prod. No. ALX-430-159) but does not inhibit the hedgehog (shh) pathway. May be used as a negative control for jervine (Prod. No. ALX-350-378).

Citations: 9

[View Online »](#)

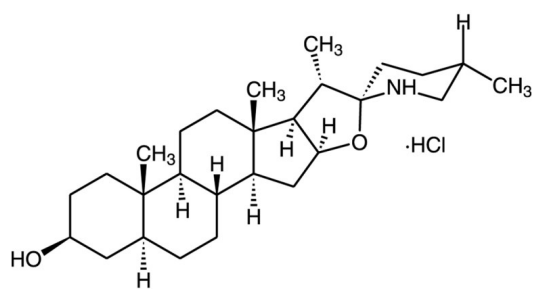
Ordering Information

[Order Online »](#)

BML-GR335-0025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C
Handling	Protect from light. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid
CAS	6192-62-7
Couple Target	Sonic hedgehog
Formula	$C_{27}H_{45}NO_2 \cdot HCl$
MW	452.1
Purity	≥90% (TLC)
Solubility	Soluble in DMSO (2mg/ml), 100% ethanol (2.5mg/ml) or dimethyl formamide (2mg/ml). Slight warming may be required.
Source	Tomato leaves

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-
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)