

Monastrol

Kinesin inhibitor

Cell permeable non-tubulin-interacting mitosis inhibitor. Blocks mitosis ($IC_{50}=14\mu M$) by binding to the mitotic kinesin Eg5.

Citations: 13

[View Online »](#)

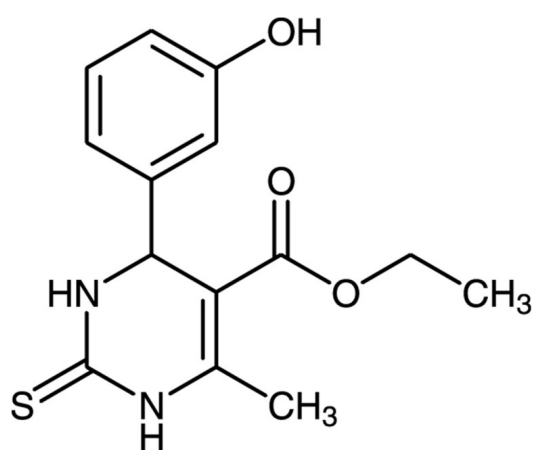
Ordering Information

[Order Online »](#)

BML-GR322-0005	5mg
BML-GR322-0025	25mg

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 -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-(3-Hydroxyphenyl)-6-methyl-2-thioxo-1,2,3,4-tetrahydro-4H-pyrimidin-5-carboxylic acid ethyl ester
Appearance	White to light yellow solid.
CAS	329689-23-8
Couple Target	Kinesin
Couple Type	Inhibitor
Formula	$C_{14}H_{16}N_2O_3S$
MW	292.4
MeltingPoint	184-186°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20mg/ml) or 100% ethanol (4mg/ml).



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)