# **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 »

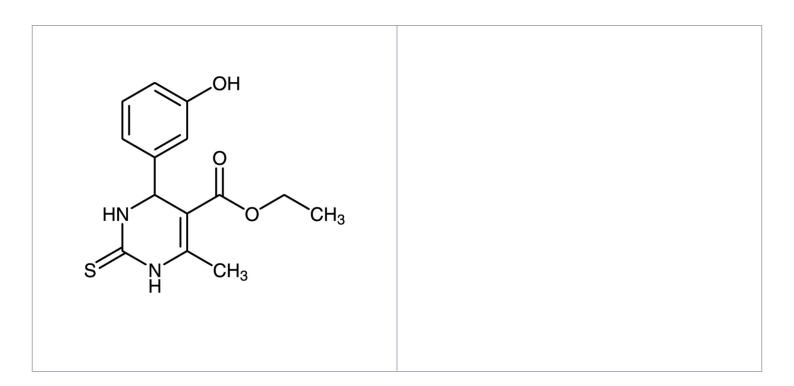
# **Ordering Information**

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)

**Soluble** in DMSO (20mg/ml) or 100% ethanol (4mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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