

ZM 336372

Raf inhibitor

Potent and specific inhibitor of the protein kinase c-Raf (IC_{50} =70nM). Has no significant effect on many other protein kinases tested (even at 50 μ M) with the exception of p38 (SAPK2a) (IC_{50} =2 μ M) and SAPK2b (p38 β) (IC_{50} =2 μ M). Induces a paradoxical >100-fold activation of c-Raf in whole cells.

Citations: 2

[View Online »](#)

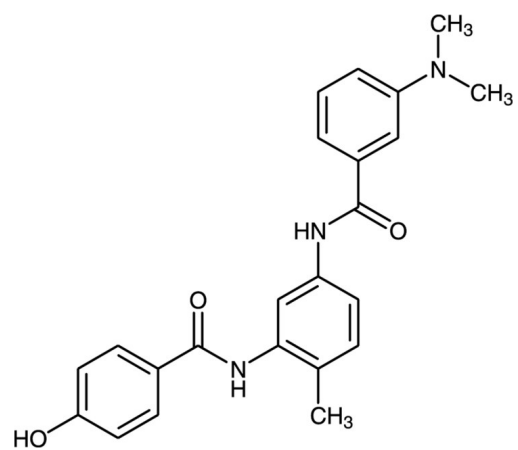
Ordering Information

[Order Online »](#)

BML-EI298-0005	5mg
BML-EI298-0025	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 3 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-[5-(3-Dimethylaminobenzamido)-2-methylphenyl]-4-hydroxybenzamide
Appearance	White to off-white solid.
CAS	208260-29-1
Couple Target	p38, Raf
Couple Type	Inhibitor
Formula	$C_{23}H_{23}N_3O_3$
MW	389.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (200mg/ml).

Technical Info / Product Notes Note: Product is not sterile.



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-be@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