

# 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 $\mu$ M) with the exception of p38 (SAPK2a) ( $IC_{50}$ =2 $\mu$ M) and SAPK2b (p38 $\beta$ ) ( $IC_{50}$ =2 $\mu$ M). Induces a paradoxical >100-fold activation of c-Raf in whole cells.

Citations: 2

[View Online »](#)

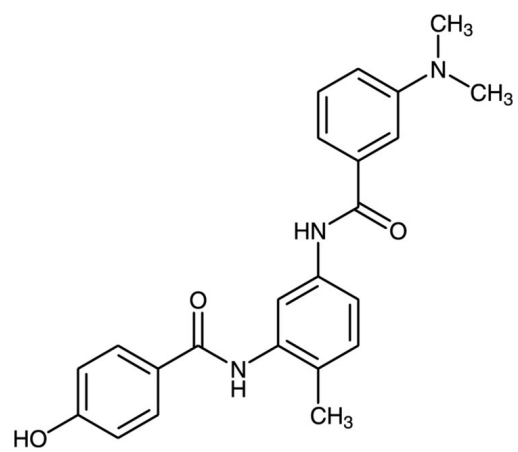
## Ordering Information

[Order Online »](#)

BML-EI298-0005	5mg
BML-EI298-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Protect from light. Keep under inert gas.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	N-[5-(3-Dimethylaminobenzamido)-2-methylphenyl]-4-hydroxybenzamide
<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	208260-29-1
<b>Couple Target</b>	p38, Raf
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{23}H_{23}N_3O_3$
<b>MW</b>	389.4
<b>Purity</b>	≥95% (HPLC)
<b>Solubility</b>	Soluble in DMSO (200mg/ml).
<b>Technical Info / Product Notes</b>	Note: Product is not sterile.

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-  
be@enzolifesciences.com](mailto:info-be@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)