

# GW5074

## Raf inhibitor

Potent and selective cell permeable inhibitor of cRAF1 kinase ( $IC_{50} = 9$  nM) with 100-fold selectivity over CDK1, CDK2, c-src, ERK2, MEK, p38, Tie2, VEGFR2 and c-fm. A useful tool for assessing the involvement of the Ras/Raf-1/ERK pathway in various signaling pathways. Displays neuroprotective effects *in vivo* through a MEK-ERK and Akt-independent mechanism.

Citations: 6

[View Online »](#)

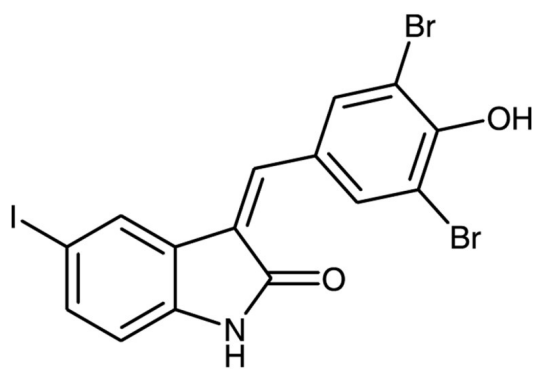
## Ordering Information

[Order Online »](#)

BML-EI307-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for at least 6 months when stored 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>	5-Iodo-3-[(3,5-dibromo-4-hydroxyphenyl)methylene]-2-indolinone
<b>Appearance</b>	Yellow to orange solid.
<b>CAS</b>	220904-83-6
<b>Couple Target</b>	Raf
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{15}H_8Br_2INO_2$
<b>MW</b>	520.9
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (>25 mg/ml).

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)