

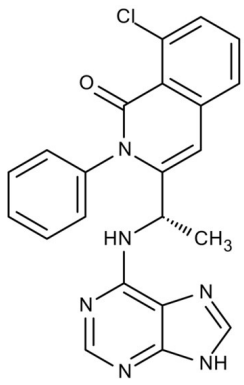
# Duvelisib (free base)

Highly selective inhibitor of PI3K- $\delta$  and PI3K- $\gamma$

Duvelisib is an orally bioavailable and highly selective inhibitor of PI3K- $\delta$  and PI3K- $\gamma$  with  $IC_{50}$  values of 0.36 nM and 19.6 nM, respectively.

**Manuals, SDS & CofA**

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	White solid.
<b>CAS</b>	1201438-56-3
<b>Couple Target</b>	PI3 kinase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{22}H_{17}ClN_6O$
<b>MW</b>	416.9
<b>Purity</b>	≥99% (HPLC)
<b>Solubility</b>	Soluble in DMSO.

**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)