

# AS1842856

## Foxo1 inhibitor

AS1842856 is a cell-permeable compound that specifically and potently inhibits the transcriptional activity of the Forkhead box O1 (Foxo1) protein (IC<sub>50</sub> value of 33 nM). This inhibition is mediated by the direct binding of AS1842856 to the active form of Foxo1, but not to its Ser256-phosphorylated (inactive) form.

AS1842856 has been shown to preferentially inhibit Foxo1 over other related proteins such as Foxo3a and Foxo4. This selectivity makes it applicable for investigating the role of Foxo1 in various cellular processes, like metabolism, cell cycle regulation, and apoptosis. Moreover, AS1842856 can inhibit gluconeogenesis in liver cells and reduce glucose production in both in vitro and in vivo models, extending its use in the research of metabolic disorders, such as diabetes.

## Ordering Information

[Order Online »](#)

ENZ-CHM325-0050	50mg
ENZ-CHM325-0100	100mg

## Manuals, SDS & CofA

[View Online »](#)

- GMP format available

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 6 months or -20°C for 1 month.
<b>Handling</b>	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	FOXO1 Inhibitor
<b>Appearance</b>	White to light yellow solid.
<b>CAS</b>	836620-48-5
<b>Couple Target</b>	FOX
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{18}H_{22}FN_3O_3$
<b>MW</b>	347.38
<b>Purity</b>	≥99% (HPLC)
<b>Solubility</b>	Soluble in DMSO (5 mg/ml).



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