

WS6

Stimulates pancreatic β -cell proliferation

WS6 is a small-molecule inhibitor of the IKB kinase pathway and ErbB3-binding protein 1 (EBP1), also known as proliferation-associated protein 2G4 (PA2G4), an RNA-binding protein involved in growth regulation.

Key features and applications include:

- **Diabetes Research:** WS6 is primarily used to study β cell proliferation, which is crucial for developing treatments for diabetes. It has been shown to increase β cell mass and normalize blood glucose levels in diabetic mouse models.
- **Endocrinology and Metabolism:** Researchers use WS6 to explore its effects on both α and β cell proliferation in human pancreatic islets, making it valuable for studies on endocrine function and metabolic diseases.
- **Cell Proliferation Studies:** WS6 is employed to investigate the mechanisms of cell proliferation, particularly through the IKB kinase pathway and ErbB3-binding protein 1 (EBP1).

Relevant disease states include:

- **Type 1 Diabetes:** WS6's ability to promote β cell growth makes it relevant for research into potential therapies for Type 1 diabetes, where β cell loss is a primary concern.
- **Metabolic Disorders:** Its role in enhancing pancreatic cell proliferation is also significant for studying various metabolic disorders that involve pancreatic dysfunction.

Ordering Information

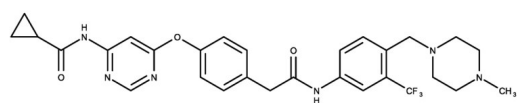
[Order Online »](#)

ENZ-CHM340-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 3 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-[[[6-[Cyclopropylcarbonyl]amino]-4-pyrimidinyl]oxy]-N-[4-[(4-methyl-1-piperazinyl)methyl]-3-(trifluoromethyl)phenyl]-benzeneacetamide
Appearance	Off-white to beige solid.
CAS	1421227-53-3
Couple Target	ErbB3-binding protein 1, IKK
Couple Type	Inhibitor
Formula	$C_{29}H_{31}F_3N_6O_3$
Identity	Determined by NMR.
MW	568.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (> 30 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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)