

# Glucagon receptor antagonist I

Potent and selective cell-permeable competitive antagonist of the glucagon receptor.

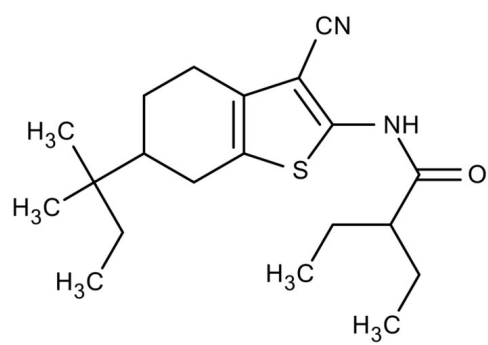
## Ordering Information

[Order Online »](#)

|                 |     |
|-----------------|-----|
| ENZ-CHM127-0005 | 5mg |
|-----------------|-----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                    |  |
|--------------------|--|
| Use/Stability      | As indicated on product label or CoA when stored as recommended. |
| Short Term Storage | -20°C  |
| Long Term Storage  | -80°C  |
| Shipping           | Ambient Temperature  |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                  |   |
|------------------|---|
| Alternative Name | hGCGR antagonist, N-[3-cyano-6-(2-methylbutan-2-yl)-4,5,6,7-tetrahydro-1-benzothiophen-2-yl]-2-ethylbutanamide, N-(3-Cyano-6-(1,1-dimethylpropyl)-4,5,6,7-tetrahydro-1-benzothien-2-yl)-2-ethylbutanamide |
| Appearance       | White crystalline solid.  |
| CAS              | 438618-32-7   |
| Couple Target    | Glucagon receptor   |
| Couple Type      | Inhibitor   |
| Formula          | $C_{20}H_{30}N_2OS$   |
| MW               | 346.5   |
| Purity           | ≥99% (HPLC)   |
| Solubility       | Soluble in DMSO or methanol.  |

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)