

# PSB-1584

## GPR84 agonist

GPR84 is an orphan G protein-coupled receptor that is activated by medium-chain fatty acids and hydroxyfatty acids although these may not be its endogenous agonist. It is highly expressed on immune cells and expression levels rise after exposure to inflammatory stimuli. Potent synthetic agonists enhance proinflammatory signaling and macrophage effector functions. PSB-1584 is a potent synthetic exogenous agonist at GPR84, which is selective against other free fatty acid receptors and is metabolically stable when incubated with human liver microsomes. An important new tool for studying the physiology of GPR84.

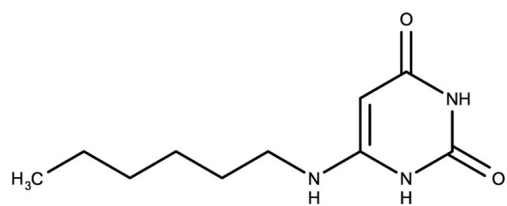
### Ordering Information

[Order Online »](#)

|                 |      |
|-----------------|------|
| ENZ-CHM479-0025 | 25mg |
|-----------------|------|

### 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** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 6-hexylamino-2,4(1H,3H)-pyrimidinedione

**Appearance** Off-white solid.

**CAS** 72255-76-6

**Couple Target** GPR84

**Couple Type** Activator

**Formula**  $C_{10}H_{17}N_3O_2$

**Identity** Determined by NMR.

**MW** 211.26

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (35 mg/ml).

**Last modified: February 27, 2025**



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)