

# CL-316243 . disodium salt

**β3-adrenergic agonist / antiobesity**

Potent and selective β-3 adrenergic receptor agonist. Displays anti-obesity and anti-diabetic effects in mice. It increases brown fat thermogenesis and metabolic rate and decreases blood insulin and glucose levels. Induces functionally active mitochondrial uncoupling protein (UCP) in white fat.

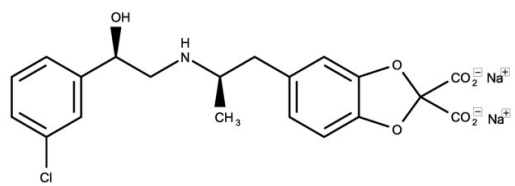
## Ordering Information

[Order Online »](#)

|                 |      |
|-----------------|------|
| ENZ-CHM483-0025 | 25mg |
|-----------------|------|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** Ambient

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5-[(2R)-2-[[[(2R)-2-(3-Chlorophenyl)-2-hydroxyethyl]amino]propyl]-1,3-benzodioxole-2,2-dicarboxylic acid disodium salt

**Appearance** Off-white solid.

**CAS** 151126-84-0

**Couple Target** Adrenergic receptor

**Couple Type** Activator

**Formula**  $C_{20}H_{18}ClNa_2O_7$

**Identity** Determined by NMR.

**MW** 465.80

**Purity**  $\geq 98\%$  (HPLC)

**Solubility** Soluble in water (50 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)