

L-755,507

β3 adrenergic agonist / CRISPR efficiency enhancer

L-755,507 is a very potent human β3 adrenergic receptor agonist. L-755,507 has been shown to increase homology-directed repair efficiency by 2- and 3-fold in CRISPR gene editing.

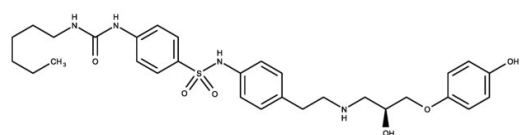
Ordering Information

[Order Online »](#)

ENZ-CHM451-0025	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 1-Hexyl-3-[4-[[4-[2-[[[(2S)-2-hydroxy-3-(4-hydroxyphenoxy)propyl]amino]ethyl]phenyl]sulfamoyl]phenyl]urea

Appearance Off-white solid.

CAS 159182-43-1

Couple Type Ligand

Formula $C_{30}H_{40}N_4O_6S$

Identity Determined by NMR.

MW 584.73

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (>25 mg/ml).

Last modified: January 24, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com