

# R-(+)-Baclofen . hydrochloride

GABA<sub>B</sub> receptor agonist

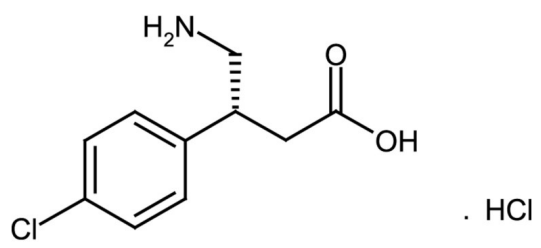
Selective GABA<sub>B</sub> receptor agonist. Skeletal muscle relaxant. The more active enantiomer of (±)-Baclofen (Prod. No. BML-EA128).

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** R-(+)-β-(Aminomethyl)-4-chlorobenzenepropanoic acid . HCl, R-(+)-4-Amino-3-(4-chlorophenyl)butanoic acid . HCl

**Appearance** White to off-white solid.

**CAS** 63701-55-3

**Couple Target** GABA receptor

**Couple Type** Activator, Ligand

**Formula**  $C_{10}H_{12}ClNO_2 \cdot HCl$

**MI** **14: 937**

**MW** 213.7 . 36.5

**Purity** ≥98% (HPLC)

**RTECS** MW5084350

**Solubility** Soluble in water, dilute aqueous base or DMSO.

**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-  
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)