

# Sertraline . HCl

Serotonin transporter inhibitor

A selective serotonin transporter inhibitor. In clinical use as an antidepressant. Displays antiproliferative activity on human colorectal carcinoma cells.

Citations: 1

[View Online »](#)

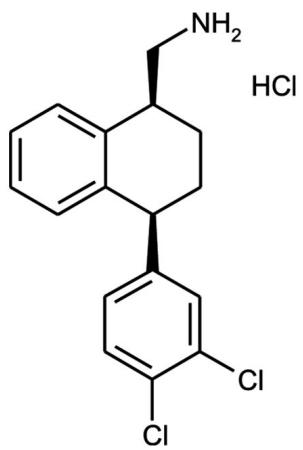
Ordering Information

[Order Online »](#)

BML-NS115-0010	10mg
----------------	------

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** CP-51974

**Appearance** White solid.

**CAS** 79559-97-0

**Couple Target** Serotonin transporter

**Couple Type** Inhibitor

**Formula**  $C_{17}H_{17}Cl_2N \cdot HCl$

**MW** 342.7

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

**Solubility** Soluble in DMSO (>25mg/ml), ethanol (10mg/ml) or water (3mg/ml).

**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)