

# Bopindolol malonate

Adrenergic receptor ligand

A slowly dissociating  $\beta_1$ -adrenergic receptor antagonist. Antihypertensive. Antioxidant ( $IC_{50}=1.8 \mu\text{M}$ ) and free radical scavenging activity.

Citations: 4

[View Online »](#)

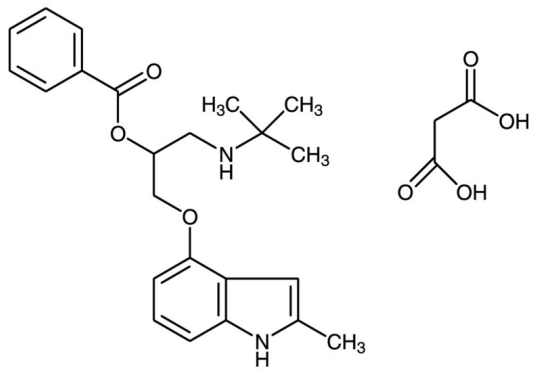
Ordering Information

[Order Online »](#)

BML-AR100-0100	100mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 4 months
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	White solid.
<b>CAS</b>	82857-38-3
<b>Couple Target</b>	Adrenergic receptor
<b>Couple Type</b>	Ligand
<b>Formula</b>	$C_{23}H_{28}N_2O_3 \cdot C_3H_4O_4$
<b>MW</b>	484.6
<b>Purity</b>	≥97% (HPLC)
<b>Solubility</b>	Soluble in water at 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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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