

Guanfacine . HCl

Adrenergic receptor ligand

Centrally acting α_2 -adrenergic receptor agonist. In the guinea pig cerebral cortex, the K_d is 20 and 880 nM for the α_{2A} - and the α_{2C} -adrenoceptor, respectively. Antihypertensive agent.

Citations: 2

[View Online »](#)

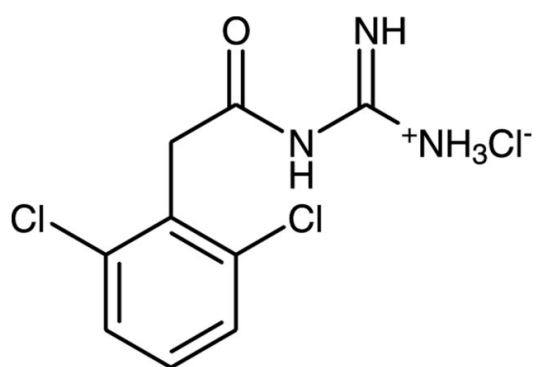
Ordering Information

[Order Online »](#)

BML-AR102-0100	100mg
BML-AR102-0500	500mg

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

Appearance White solid.

CAS 29520-14-7 / 29110-48-3

Couple Target Adrenergic receptor

Couple Type Ligand

Formula $C_9H_9Cl_2N_3O \cdot HCl$

MW 282.5

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in water (25 mg/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)