

(+/-)Cyanopindolol hemifumarate

Adrenergic receptor and serotonin receptor
antagonist

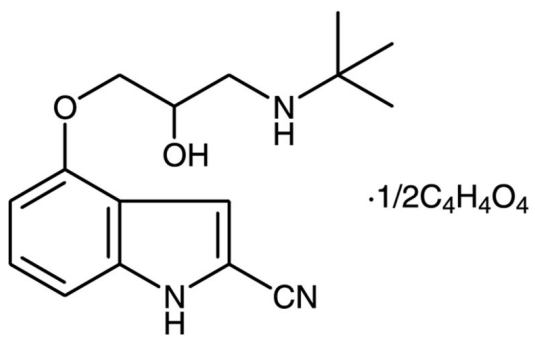
β -Adrenergic receptor antagonist. Serotonin 5HT1A/1B antagonist.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	69906-85-0
Couple Target	Adrenergic receptor, Serotonin receptor
Couple Type	Ligand
Formula	$C_{16}H_{21}N_3O_2 \cdot 1/2C_4H_4O_4$
MW	345.4
Purity	≥99% (TLC)
Solubility	Soluble in warm water (2 mg/ml), DMSO (8 mg/ml).

Last modified: May 29, 2024



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-be@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