

R-(+)-DIOA

K⁺ channel blocker

Inhibitor of the K⁺/Cl⁻-cotransport that does not affect the bumetanide-sensitive Na⁺/K⁺/Cl⁻-cotransport system.

Citations: 3

[View Online »](#)

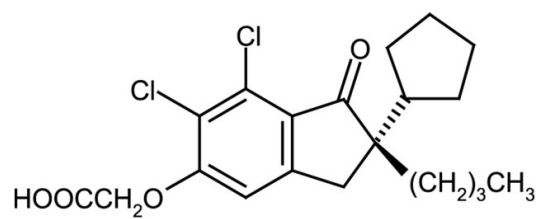
Ordering Information

[Order Online »](#)

ALX-550-263-M005	5mg
ALX-550-263-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name R-(+)-Butylindazone, (Dihydroindenyl)oxy alkanoic acid

Appearance White solid.

CAS 81166-47-4

Couple Target Potassium channel

Couple Type Blocker

Formula $C_{20}H_{24}Cl_2O_4$

MW 399.3

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol or dilute aqueous base; insoluble in water or dilute aqueous acid.

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-fr@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