

R-(-)-Glucoraphanin . potassium salt

Cytochrome P450 modulator

Modulator of hepatic cytochrome P450 and phase II conjugation activity.
The enzyme myrosinase transforms R-(-)-glucoraphanin into R-Sulforaphane (Prod. No. ALX-350-230).

Citations: 4

[View Online »](#)

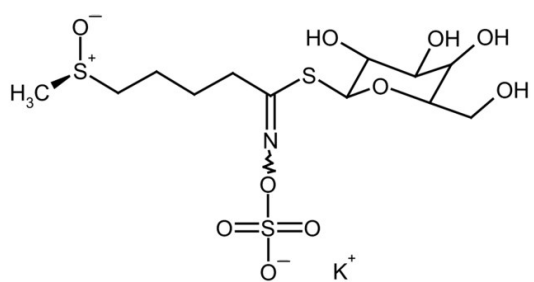
Ordering Information

[Order Online »](#)

ALX-350-181-M010	10mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Very hygroscopic.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-Methylsulfinylbutylglucosinate . K
Appearance	White solid.
CAS	21414-41-5
Couple Target	Cytochrome P450
Formula	$C_{12}H_{22}NO_{10}S_3 \cdot K \cdot xH_2O$
MW	436.5 . 39.1 (anhydrous basis)
Purity	≥99% (HPLC)
Source	Isolated from <i>Brassica oleracea</i> var. <i>palmifolia</i>

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-
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)