

Ruthenium red

Capsaicin and Ca^{2+} antagonist

Blocks TRPV1. Capsaicin and calcium antagonist. Inhibitor of $\text{Ca}^{2+}/\text{Mg}^{2+}$ -ATPase.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-550-295-G001	1g
------------------	----

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 Dark red to purple crystalline solid.

CAS 11103-72-3

Formula $\text{Cl}_6\text{H}_{42}\text{N}_{14}\text{O}_2\text{Ru}_3 \cdot 4\text{H}_2\text{O}$

MI **14:** 8300

MW 786.4 . 72.1

Purity $\geq 90\%$ (metals basis)

RTECS VM2656500

Sequence $[(\text{NH}_3)_5\text{RuORu}(\text{NH}_3)_4\text{ORu}(\text{NH}_3)_5]^{6+} 6\text{Cl}^-$

Solubility Soluble in water (10mg/ml) or ammonia.

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)