

RITA

Inhibitor of p53/HDM2 interaction

Inhibitor of p53/HDM2 interaction. Activator of p53 function in tumors. Potently inhibits the growth of a variety of cancer cell lines in the submicromolar range. It has been suggested that RITA selectively targets cells expressing p53 by other mechanisms than direct binding to the N-terminal domain of p53.

Citations: 4

[View Online »](#)

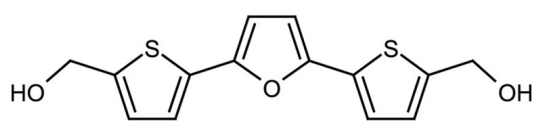
Ordering Information

[Order Online »](#)

ALX-430-151-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least two years after receipt when stored at -20°C. We do not recommend storing aqueous solutions for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Reactivation of p53 and Induction of Tumor Cell Apoptosis, 2,5-bis(5-Hydroxymethyl-2-thienyl)furan, NSC652287
Appearance	White to brown crystalline solid.
CAS	213261-59-7
Formula	$C_{14}H_{12}O_3S_2$
MW	292.4
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol (1mg/ml), DMSO (20mg/ml), dimethyl formamide (30mg/ml); 0.5mg/ml soluble in a 1:1 solution of dimethyl formamide: PBS (pH 7.2) (dissolve RITA first in dimethyl formamide and then dilute with PBS).



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)