

t-Butylhydroquinone

Ca²⁺ ATPase inhibitor. Erk2 activator.

Inhibitor of the endoplasmic reticulum Ca²⁺-ATPase. Activates extracellular signal-regulated protein kinase 2 (ERK2).

Citations: 1

[View Online »](#)

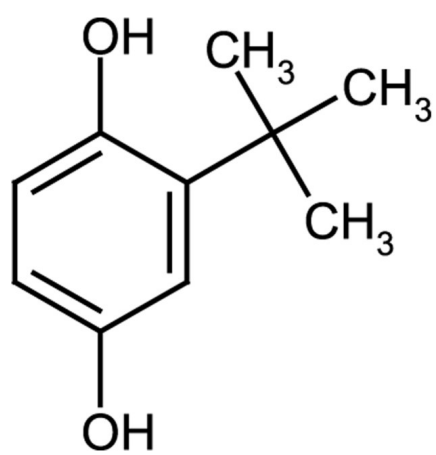
Ordering Information

[Order Online »](#)

ALX-270-061-G010	10g
------------------	-----

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

Alternative Name tBHQ

Appearance White to off-white solid.

CAS 1948-33-0

Couple Target ATPase, Erk

Couple Type Activator, Inhibitor

Formula $C_{10}H_{14}O_2$

Identity Determined by IR.

MW 166.2

Purity $\geq 97\%$ (GC)

Solubility Soluble in DMSO or methanol.

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)