

Hydroxytyrosol . acetate

Antioxidant

Antioxidant comparable to resveratrol (Prod. No. BML-FR104). Is converted to hydroxytyrosol through hydrolysis by esterases. Reaches full oxygen radical absorbance capacity (ORAC) when converted to free hydroxytyrosol. ORAC values of hydroxytyrosol are about 50% higher than resveratrol, 10-fold higher than L(+)-ascorbic acid and 4-fold higher than (-)-epicatechin. Has been shown to have anticancer and antiviral effects. Antibacterial at 400µg/mL against *Streptococcus pyogenes*, *Staphylococcus aureus*, *Escherichia coli* and *Klebsiella pneumoniae*.

Hydroxytyrosol-acetate is a small-molecule phenolic compound found in the leaves and fruits of olive (*Olea europaea*). It is a metabolite of oleuropein (Ole), one of the major polyphenolic components of olive products.

Citations: 8

[View Online »](#)

Ordering Information

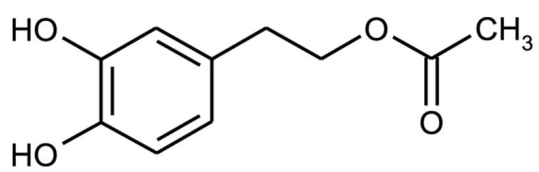
[Order Online »](#)

ALX-350-404-M050

50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HTS, 2-(3,4-Dihydroxyphenyl)-ethyl acetate
Appearance	White to off-white powder.
CAS	69039-02-7
Formula	$C_{10}H_{12}O_4$
MW	196.2
Purity	≥98% (HPLC)
Solubility	Freely soluble in organic solvents; slightly soluble in water (10mg/ml).
Source	Synthetic

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)