

# Hispolon

## Apoptosis inducer

Polyphenolic compound. Induces ROS-mediated apoptosis. Has antiviral activity.

Citations: 5

[View Online »](#)

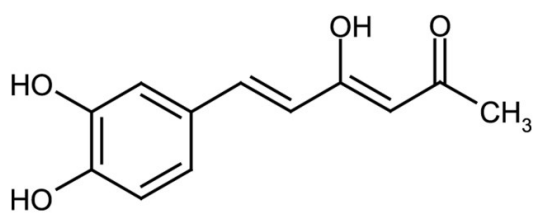
### Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-350-384-M005 | 5mg  |
| ALX-350-384-M025 | 25mg |

### 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

|            |                    |
|------------|--------------------|
| Appearance | Yellow solid.      |
| CAS        | 173933-40-9        |
| Formula    | $C_{12}H_{12}O_4$  |
| Identity   | Determined by NMR. |
| MW         | 220.2              |
| Purity     | ≥98% (HPLC)        |
| Solubility | Soluble in DMSO.   |
| 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-  
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)