

# Cholesterol 3-sulfate . sodium salt

**PKC activator**

Activates multiple protein kinase C (PKC) iso-enzymes, especially the  $\epsilon$ ,  $\eta$  and  $\zeta$  isoforms.

Citations: 2

[View Online »](#)

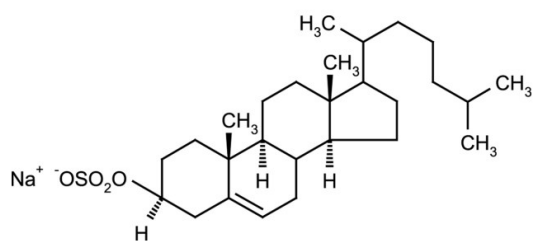
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-370-009-M005 | 5mg |
|------------------|-----|

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 | +4°C                                                             |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|               |                             |
|---------------|-----------------------------|
| Appearance    | White to off-white powder.  |
| CAS           | 2864-50-8                   |
| Couple Target | PKC                         |
| Couple Type   | Activator                   |
| Formula       | $C_{27}H_{45}O_4S \cdot Na$ |
| Identity      | Determined by IR.           |
| MW            | 465.7 . 23.0                |
| Purity        | One spot by TLC.            |
| Solubility    | Soluble in 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)