

# 6-ECDCA

## FXR ligand

A novel derivative of cholic acid which acts as a potent and selective FXR agonist ( $EC_{50} = 99\text{nM}$ ) displaying anticholeretic activity in an *in vivo* rat model of cholestasis. It not only inhibits vascular smooth muscle cell inflammation and migration but also promotes adipocyte differentiation and regulates adipose cell function *in vivo*.

Citations: 4

[View Online »](#)

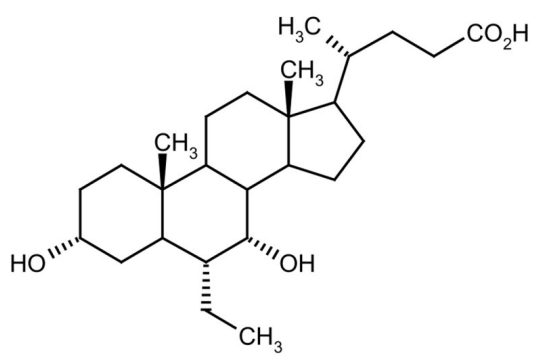
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-GR246-0005 | 5mg  |
| BML-GR246-0025 | 25mg |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |   |
|-------------------|---|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months. |
| Long Term Storage | -20°C   |
| Shipping          | Blue Ice  |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |   |
|------------------|---|
| Alternative Name | 6-Ethylchenodeoxycholic acid, 3 $\alpha$ ,7 $\alpha$ -Dihydroxy-6 $\alpha$ -ethyl-5 $\beta$ -cholan-24-oic acid, Obeticholic acid |
| Appearance       | White solid.  |
| CAS              | 459789-99-2   |
| Couple Target    | FXR   |
| Couple Type      | Ligand  |
| Formula          | C <sub>26</sub> H <sub>44</sub> O <sub>4</sub>  |
| MW               | 420.6   |
| Purity           | ≥97% (TLC)  |
| Solubility       | Soluble in DMSO (35mg/ml) or 100% ethanol (25mg/ml).  |

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)