

# Forskolin

## Adenylate cyclase activator

Forskolin is naturally produced by the Coleus plant, *Coleus forskohlii*. Forskolin activates adenylate cyclase by directly interacting with the catalytic unit of the enzyme leading to an increase in the intracellular concentration of cAMP. Forskolin is a widely used tool for the investigation of the role of cAMP as a second messenger with a broad range of potential therapeutic applications. Forskolin is a ionotropic agent and vasodilator. Forskolin also reduces platelet aggregation in a dose-dependent manner, inhibits ion channels by a mechanism that does not involve cAMP, inhibits nicotinic acetylcholine receptors and induces neuronal differentiation in stem cells and several neuroblastoma.

Citations: 45

[View Online »](#)

## Ordering Information

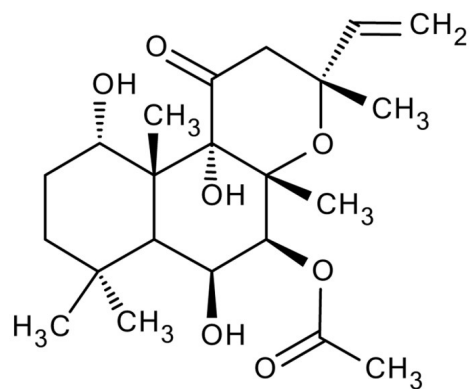
[Order Online »](#)

|                |       |
|----------------|-------|
| BML-CN100-0010 | 10mg  |
| BML-CN100-0100 | 100mg |

## Manuals, SDS & CofA

[View Online »](#)

- Cell-permeable, potent, reversile, and rapid activator of adenylate cyclase
- Compound used to study the role of cAMP as secondary messenger
- Highly cited



## Handling & Storage

|                   |   |
|-------------------|---|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for up to 4 months when stored at +4°C. Do not dissolve in ethanol, which inhibits forskolin activation of adenylate cyclase. |
| Long Term Storage | -20°C   |
| Shipping          | Ambient Temperature   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |  |
|------------------|--|
| Alternative Name | Coleonol, Colforsin  |
| Appearance       | White solid.   |
| CAS              | 66575-29-9   |
| Couple Target    | Adenylate cyclase  |
| Couple Type      | Activator  |
| Formula          | $C_{22}H_{34}O_7$  |
| MI               | 14: 2476   |
| MW               | 410.5  |
| Purity           | ≥97% (HPLC, TLC)   |
| Solubility       | Soluble in anhydrous DMSO (5mg/ml), 100% ethanol (6mg/ml), or ethyl acetate (10mg/ml). DMSO is the recommended solvent for activation of cAMP. |
| Source           | Isolated from <i>Coleus forskohlii</i> .   |

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)