

# RO-20-1724

## PDE inhibitor

Selective, potent cAMP-specific phosphodiesterase inhibitor (PDE IV,  $IC_{50}$  =2 $\mu$ M). Inhibits superoxide generation and platelet aggregation stimulated by arachidonic acid. Also inhibits fMLP-induced neutrophil adhesion to vascular endothelial cells.

Citations: 10

[View Online »](#)

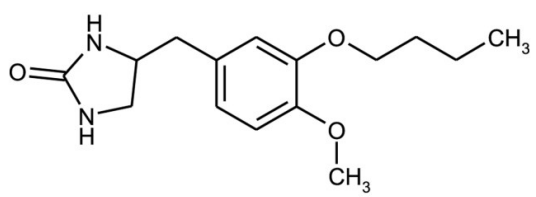
## Ordering Information

[Order Online »](#)

|                |       |
|----------------|-------|
| BML-EI117-0100 | 100mg |
| BML-EI117-1000 | 1g    |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                                                                                                                                               |
|-------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Solutions are stable for up to 3 months when stored at -20°C. |
| Long Term Storage | Ambient                                                                                                                                                                                       |
| Shipping          | Ambient Temperature                                                                                                                                                                           |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                            |
|------------------|----------------------------------------------------------------------------|
| Alternative Name | 4-(3-Butoxy-4-methoxybenzyl)-2-imidazolidinone                             |
| Appearance       | White to off-white solid.                                                  |
| CAS              | 29925-17-5                                                                 |
| Couple Target    | PDE                                                                        |
| Couple Type      | Inhibitor                                                                  |
| Formula          | $C_{15}H_{22}N_2O_3$                                                       |
| MW               | 278.3                                                                      |
| MeltingPoint     | 126-127°C                                                                  |
| Purity           | ≥99% (TLC)                                                                 |
| Solubility       | Soluble in 100% ethanol (40mg/ml) or DMSO (200 mg/ml); insoluble in water. |

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)