

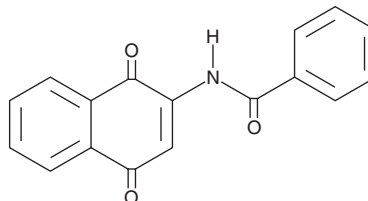
# PRODUCT INFORMATION



## PPM-18

Item No. 13327

**CAS Registry No.:** 65240-86-0  
**Formal Name:** N-(1,4-dihydro-1,4-dioxo-2-naphthalenyl)-benzamide  
**Synonym:** NSC 73233  
**MF:** C<sub>17</sub>H<sub>11</sub>NO<sub>3</sub>  
**FW:** 277.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 257, 282 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

PPM-18 is supplied as a crystalline solid. A stock solution may be made by dissolving the PPM-18 in the solvent of choice, which should be purged with an inert gas. PPM-18 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PPM-18 in these solvents is approximately 3 and 5 mg/ml, respectively.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

PPM-18 is a naphthoquinone compound that inhibits NF-κB activation *in vitro* and *in vivo* (IC<sub>50</sub> = 5 μM).<sup>1,2</sup> Through this action, it inhibits the expression of inducible nitric oxide synthase (iNOS) in macrophages treated with lipopolysaccharide and suppresses inflammation and sepsis.<sup>1</sup>

### References

1. Yu, S.M., Wu, J.F., Lin, T.L., *et al.* Inhibition of nitric oxide synthase expression by PPM-18, a novel anti-inflammatory agent, *in vitro* and *in vivo*. *Biochem. J.* **328**, 363-369 (1997).
2. Davis, M.E., Grumbach, I.M., Fukai, T., *et al.* Shear stress regulates endothelial nitric-oxide synthase promoter activity through nuclear factor κB binding. *J. Biol. Chem.* **279**(1), 163-168 (2004).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/01/2021

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)