

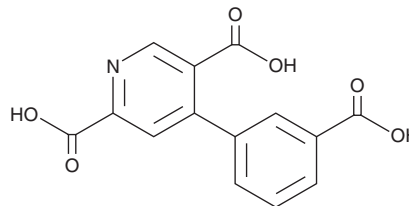
# PRODUCT INFORMATION



## 4-CPPC

Item No. 30068

**CAS Registry No.:** 29553-70-6  
**Formal Name:** 4-(3-carboxyphenyl)-2,5-pyridinedicarboxylic acid  
**MF:** C<sub>14</sub>H<sub>9</sub>NO<sub>6</sub>  
**FW:** 287.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 223 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

4-CPPC is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-CPPC in the solvent of choice, which should be purged with an inert gas. 4-CPPC is soluble in the organic solvent DMSO.

### Description

4-CPPC is an inhibitor of macrophage migration inhibitory factor-2 (MIF-2; IC<sub>50</sub> = 27 μM).<sup>1</sup> It is selective for MIF-2 over MIF-1 (IC<sub>50</sub> = 450 μM). It inhibits MIF-2, but not MIF-1, binding to CD74 *in vitro* when used at a concentration of 10 μM. 4-CPPC (5, 10, and 25 μM) inhibits MIF-2-induced ERK1/2 phosphorylation in primary human skin fibroblasts.

### Reference

1. Tilstam, P.V., Pantouris, G., Corman, M., *et al.* A selective small-molecule inhibitor of macrophage migration inhibitory factor-2 (MIF-2), a MIF cytokine superfamily member, inhibits MIF-2 biological activity. *J. Biol. Chem.* **294**(49), 18522-18531 (2019).

#### 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, 04/07/2020

#### 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