

# PRODUCT INFORMATION



**MK-0893**

Item No. 28923

**CAS Registry No.:** 870823-12-4  
**Formal Name:** N-[4-[(1S)-1-[3-(3,5-dichlorophenyl)-5-(6-methoxy-2-naphthalenyl)-1H-pyrazol-1-yl]ethyl]benzoyl]-β-alanine

**MF:** C<sub>32</sub>H<sub>27</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>4</sub>

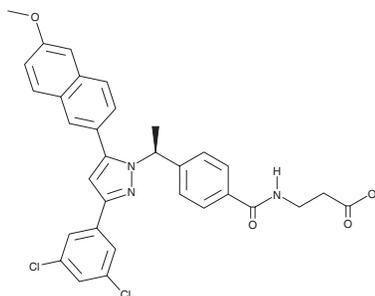
**FW:** 588.5

**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years



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

## Laboratory Procedures

MK-0893 is supplied as a solid. A stock solution may be made by dissolving the MK-0893 in the solvent of choice, which should be purged with an inert gas. MK-0893 is soluble (≥10 mg/ml) in ethanol and DMSO.

## Description

MK-0893 is an allosteric antagonist of the glucagon receptor (GCGR; IC<sub>50</sub> = 15.7 nM).<sup>1,2</sup> It is selective for GCGR over glucose-dependent insulinotropic polypeptide receptor (GIPR), glucagon-like peptide 1 receptor (GLP-1R), pituitary adenylate cyclase-activating peptide (PACAP) type 1 (PAC<sub>1</sub>) receptor, vasoactive intestinal polypeptide receptor 1 (VPAC<sub>1</sub>), and VPAC<sub>2</sub> (IC<sub>50</sub>s = 1,020, >10,000, 9,200, 11,580, and 12,800 nM, respectively).<sup>2</sup> MK-0893 (1 and 3 mg/kg) reduces glucagon-induced increases in blood glucose levels in rhesus monkeys.

## References

1. Jazayeri, A., Doré, A.S., Lamb, D., *et al.* Extra-helical binding site of a glucagon receptor antagonist. *Nature* **533(7602)**, 274-277 (2016).
2. Xiong, Y., Guo, J., Candelore, M.R., *et al.* Discovery of a novel glucagon receptor antagonist N-[(4-[(1S)-1-[3-(3,5-dichlorophenyl)-5-(6-methoxynaphthalen-2-yl)-1H-pyrazol-1-yl]ethyl]phenyl)carbonyl]-β-alanine (MK-0893) for the treatment of type II diabetes. *J. Med. Chem.* **55(13)**, 6137-6148 (2012).

### 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, 02/04/2026

## 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