

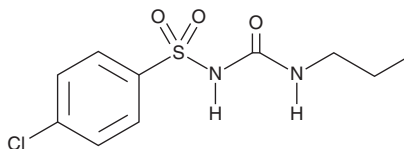
# PRODUCT INFORMATION



## Chlorpropamide

Item No. 26087

**CAS Registry No.:** 94-20-2  
**Formal Name:** 4-chloro-N-[(propylamino) carbonyl]-benzenesulfonamide  
**Synonyms:** NSC 44634, NSC 626720  
**MF:** C<sub>10</sub>H<sub>13</sub>ClN<sub>2</sub>O<sub>3</sub>S  
**FW:** 276.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232 nm  
**Supplied as:** A crystalline solid  
**Storage:** 4°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Chlorpropamide is supplied as a crystalline solid. A stock solution may be made by dissolving the chlorpropamide in the solvent of choice, which should be purged with an inert gas. Chlorpropamide is soluble in organic solvents such as ethanol, acetone, and chloroform.

### Description

Chlorpropamide is a first generation sulfonyleurea.<sup>1</sup> It increases glycolysis by increasing fructose-2,6-bisphosphate levels and inhibits glucagon-induced gluconeogenesis by augmenting the effect of insulin in isolated rat hepatocytes when used at a concentration of 0.2 mM. Chlorpropamide reverses increases in blood glucose levels induced by hydrochlorothiazide (Item No. 21304) and previous insulin in a rabbit model of diabetes when administered in a single 250 mg/kg dose or repeated doses of 10 mg/kg per day for six days.<sup>2</sup> Formulations containing chlorpropamide have been used as adjuncts to diet and exercise in the treatment of type 2 diabetes.

### References

1. Monge, L., Mojena, M., Ortega, J.L., *et al.* Chlorpropamide raises fructose-2,6-bisphosphate concentration and inhibits gluconeogenesis in isolated rat hepatocytes. *Diabetes* **35(1)**, 89-96 (1986).
2. Pakrashi, A. and Mukherjee, S.K. A new method for the production of experimental diabetes in rabbits. *J. Endocrinol.* **66(2)**, 147-150 (1975).

#### 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, 11/07/2022

#### 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