

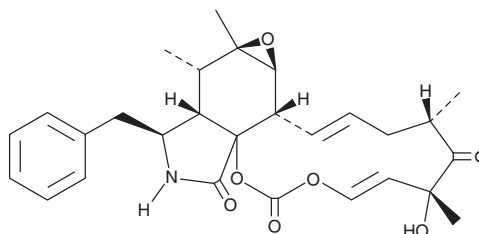
# PRODUCT INFORMATION



**Gö 6983**

Item No. 13311

**CAS Registry No.:** 133053-19-7  
**Formal Name:** 3-[1-[3-(dimethylamino)propyl]-5-methoxy-1H-indol-3-yl]-4-(1H-indol-3-yl)-1H-pyrrole-2,5-dione  
**MF:** C<sub>26</sub>H<sub>26</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 442.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

Gö 6983 is supplied as a crystalline solid. A stock solution may be made by dissolving the Gö 6983 in an organic solvent purged with an inert gas. Gö 6983 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of Gö 6983 in these solvents is approximately 10 mg/ml.

Gö 6983 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Gö 6983 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Gö 6983 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Gö 6983 inhibits several isoforms of protein kinase C (PKC; IC<sub>50</sub> = 7, 7, 6, 10, 60, and 20,000 nM for PKC $\alpha$ , PKC $\beta$ , PKC $\gamma$ , PKC $\delta$ , PKC $\zeta$ , and PKC $\mu$ , respectively).<sup>1</sup> It provides cardioprotective effects in myocardial ischemia/reperfusion by attenuating contractile dysfunction.<sup>2,3</sup>

## References

1. Gschwendt, M., Dieterich, S., Rennecke, J., *et al.* Inhibition of protein kinase C  $\mu$  by various inhibitors. Differentiation from protein kinase c isoenzymes. *FEBS Lett.* **392**, 77-80 (1996).
2. Peterman, E.E., Taormina, P.II., Harvey, M., *et al.* Gö 6983 exerts cardioprotective effects in myocardial ischemia/reperfusion. *J. Cardiovasc. Pharmacol.* **43**, 645-656 (2004).
3. Young, L.H., Balin, B.J., and Weis, M.T. Gö 6983: A fast acting protein kinase C inhibitor that attenuates myocardial ischemia/reperfusion injury. *Cardiovascular Drug Reviews* **23(3)**, 255-272 (2005).

### 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, 11/18/2022

## 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