

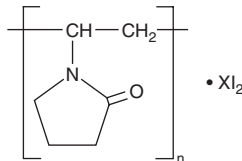
PRODUCT INFORMATION



Povidone-Iodine

Item No. 27883

CAS Registry No.: 25655-41-8
Formal Name: 1-ethenyl-2-pyrrolidinone, homopolymer, compd. with iodine
Synonyms: Iodopovidone, NSC 26245, Polyvinylpyrrolidone-Iodine
MF: $(C_6H_9NO)_n \cdot XI_2$
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Povidone-iodine is supplied as a solid. A stock solution may be made by dissolving the povidone-iodine in the solvent of choice. Povidone-iodine is soluble in the organic solvent DMSO, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Povidone-iodine is a microcide with broad-spectrum activity against bacteria, fungi, viruses, and protozoans.¹ It reduces total anaerobic and aerobic bacterial counts *in vitro* in dog feces when applied topically as a 10% solution.² Povidone-iodine decreases bacterial counts and the occurrence of anastomotic leaks in a dog model of emergent resection and anastomosis of the sigmoid colon when applied *in situ* at a concentration of 10% v/v. Topical administration of povidone-iodine (10% v/v) also reduces bacterial counts on the skin of cynomolgus monkeys.³ Formulations containing povidone-iodine have been used to disinfect skin before and after surgery.

References

1. Eggers, M., Eickmann, M., Kowalski, K., *et al.* Povidone-iodine hand wash and hand rub products demonstrated excellent *in vitro* virucidal efficacy against Ebola virus and modified vaccinia virus Ankara, the new European test virus for enveloped viruses. *BMC Infect. Dis.* **15:375** (2015).
2. Jones, F.E., DeCosse, J.J., and Condon, R.E. Evaluation of "instant" preparation of the colon with povidone-iodine. *Ann. Surg.* **184(1)**, 74-79 (1976).
3. Nakata, H., Tsubotani, Y., Nii, T., *et al.* Effects of olanexidine gluconate on preoperative skin preparation: An experimental study in cynomolgus monkeys. *J. Med. Microbiol.* **66(5)**, 678-685 (2017).

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, 12/09/2022

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