

# PRODUCT INFORMATION



## Halometasone

Item No. 34848

CAS Registry No.: 50629-82-8

Formal Name: (6 $\alpha$ ,11 $\beta$ ,16 $\alpha$ )-2-chloro-6,9-difluoro-11,17,21-trihydroxy-16-methylpregna-1,4-diene-3,20-dione

Synonyms: Ba 48401, 2-Chloroflumethasone, C 48401Ba, 2-chloro Flumethasone, Halomethasone

MF: C<sub>22</sub>H<sub>27</sub>ClF<sub>2</sub>O<sub>5</sub>

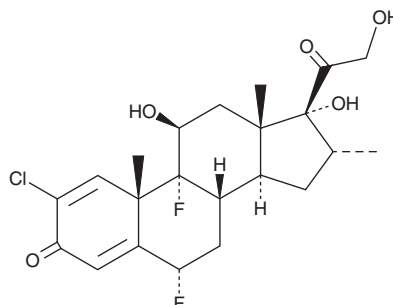
FW: 444.9

Purity:  $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Halometasone is supplied as a solid. A stock solution may be made by dissolving the halometasone in the solvent of choice, which should be purged with an inert gas. Halometasone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of halometasone in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Halometasone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, halometasone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Halometasone has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Halometasone is a synthetic corticosteroid.<sup>1,2</sup> Formulations containing halometasone have been used in the treatment of psoriasis vulgaris and eczematous dermatoses.

### References

1. de la Brassine, M., Kint, A., Lachapelle, J.M., *et al.* Halomethasone (C 48.401-Ba) for the topical treatment of common dermatoses. *J. Int. Med. Res.* **12**(5), 307-309 (1984).
2. Zhu, J.-W., Wu, X.-J., Lu, Z.-F., *et al.* Role of VEGF receptors in normal and psoriatic human keratinocytes: Evidence from irradiation with different UV sources. *PLoS One* **8**(1), e55463 (2013).

#### 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, 12/12/2022

#### 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