

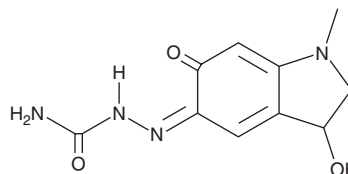
# PRODUCT INFORMATION



## Carbazochrome

Item No. 32951

**CAS Registry No.:** 69-81-8  
**Formal Name:** 2-(1,2,3,6-tetrahydro-3-hydroxy-1-methyl-6-oxo-5H-indol-5-ylidene)-hydrazinecarboxamide  
**Synonyms:** Adrenochrome Monosemicarbazone, Adrenochrome Semicarbazone, NSC 73742  
**MF:** C<sub>10</sub>H<sub>12</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 236.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 358 nm  
**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

Carbazochrome is supplied as a crystalline solid. A stock solution may be made by dissolving the carbazochrome in the solvent of choice, which should be purged with an inert gas. Carbazochrome is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of carbazochrome in these solvents is approximately 1 mg/ml.

Carbazochrome is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, carbazochrome should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Carbazochrome has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Carbazochrome is a hemostatic agent.<sup>1,2</sup> *In vivo*, carbazochrome (0.1 mg/kg) reduces blood loss by 45% in a mouse model of traumatic hemorrhage.<sup>1</sup> Topical administration of carbazochrome induces blood coagulation at back wound sites in mice.<sup>2</sup>

### References

1. Keith, J.H., Keith, W.W., and Nelson, M.H. Effects of adrenochrome semicarbazone on blood loss in the mouse. *J. Pharm. Sci.* **56**(8), 997-1001 (1967).
2. Mukherjee, D. and Banthia, A.K. Preparation of adrenochrome hydrogel patch, gel, ointment, and the comparison of their blood coagulating and wound healing capability. *Mater. Manuf. Process.* **21**(3), 297-301 (2006).

#### 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