

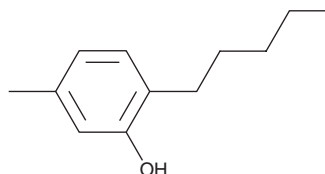
# PRODUCT INFORMATION



## Amylmetacresol

Item No. 40290

CAS Registry No.: 1300-94-3  
Formal Name: 5-methyl-2-pentyl-phenol  
Synonyms: AMC, 6-*n*-Amyl-*m*-cresol  
MF: C<sub>12</sub>H<sub>18</sub>O  
FW: 178.3  
Purity: ≥98%  
Supplied as: A neat liquid  
Storage: -20°C  
Stability: ≥4 years



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

### Laboratory Procedures

Amylmetacresol is supplied as a neat liquid. A stock solution may be made by dissolving the amylmetacresol in the solvent of choice, which should be purged with an inert gas. Amylmetacresol is soluble (≥10 mg/ml) in ethanol and DMSO.

### Description

Amylmetacresol is a topical antiseptic and analgesic agent.<sup>1</sup> It inhibits depolarization-induced whole-cell sodium currents in HEK293 cells expressing human voltage-gated sodium channel 1.2 (Na<sub>v</sub>1.2; EC<sub>50</sub> = 56.3 μM). Formulations containing amylmetacresol have been used in oral lozenges for the treatment of pharyngitis.

### Reference

1. Buchholz, V., Leuwer, M., Ahrens, J., *et al.* Topical antiseptics for the treatment of sore throat block voltage-gated neuronal sodium channels in a local anaesthetic-like manner. *Naunyn. Schmiedeberg's. Arch. Pharmacol.* **380**(2), 161-168 (2009).

#### 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, 04/05/2024

#### 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](http://WWW.CAYMANCHEM.COM)