CERTIFICATE of ANALYSIS



Methallylescaline (hydrochloride)

≥95%

Item No. 19808 * Batch No. 0810103

Purity Specification:

Molecular formula: C14H21NO3 • HCl

CAS number: 21JUL2029 Formula weight: 287.80 **Expiration date:**

Overview

Tests	Results
16513	Results
HPLC	Purity: 97.1 %
IR	Conforms
Mass Spec	M+H: 252.3
	M+ACN+H: 293.3
	2M+H: 503.4
NMR	Conforms
Counterion	Conforms
Counterior	Comornia

Reviewed and approved by: Ryan Seiffert

WARNING
THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER
TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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