

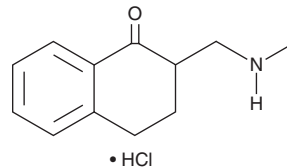
# PRODUCT INFORMATION



## MTTA (hydrochloride)

Item No. 17121

**CAS Registry No.:** 2749302-69-8  
**Formal Name:** 3,4-dihydro-2-[(methylamino)methyl]-1(2H)-naphthalenone, monohydrochloride  
**Synonym:** Mephtetramine  
**MF:** C<sub>12</sub>H<sub>15</sub>NO • HCl  
**FW:** 225.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251, 294 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

MTTA is a designer drug that features a carbonyl group adjacent to a phenyl group, similar to cathinones.<sup>1</sup> Also marketed as mephtetramine, MTTA differs structurally from cathinones, most notably by the presence of an additional carbon between the carbonyl and amine. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

### Reference

1. Power, J.D., Scott, K.R., Gardner, E.A., *et al.* The syntheses, characterization and *in vitro* metabolism of nitracaine, methoxypiperamide and mephtetramine. *Drug Test Anal.* **6(7-8)**, 668-675 (2014).

**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, 07/05/2023

### 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