

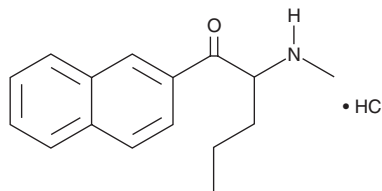
# PRODUCT INFORMATION



## NRG-3 (hydrochloride)

Item No. 14115

**CAS Registry No.:** 2701928-05-2  
**Formal Name:** 2-(methylamino)-1-(naphthalen-2-yl)pentan-1-one, monohydrochloride  
**MF:** C<sub>16</sub>H<sub>19</sub>NO • HCl  
**FW:** 277.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 208, 258, 294, 341 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.

### Laboratory Procedures

NRG-3 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the NRG-3 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. NRG-3 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NRG-3 (hydrochloride) in DMSO is approximately 2 mg/ml and approximately 1 mg/ml in ethanol and DMF.

NRG-3 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NRG-3 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NRG-3 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Substituted cathinones are psychostimulants which have appeared in chemical mixtures sold over the counter.<sup>1-3</sup> NRG-3 is a substituted cathinone that is structurally analogous to pentylone (Item No. 9000746), with the methylenedioxy group replaced with a naphthyl group. The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research applications.

### References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-15 (2011).
2. Brandt, S.D., Sumnall, H.R., Measham, F., *et al.* Analyses of second-generation 'legal-highs' in the UK: Initial findings. *Drug Test. Anal.* **2(8)**, 377-382 (2010).
3. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. Brandt, S.D., *et al.* Retrieved August 18, 2010, from [www.drugtestinganalysis.com](http://www.drugtestinganalysis.com).

#### 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, 07/14/2023

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