

# PRODUCT INFORMATION



## DOTA-NOC (acetate)

Item No. 41975

**Formal Name:** N-[2-[4,7,10-tris(carboxymethyl)-1,4,7,10-tetraazacyclododec-1-yl]acetyl]-D-phenylalanyl-L-cysteinyl-3-(1-naphthalenyl)-L-alanyl-D-tryptophyl-L-lysyl-L-threonyl-N-[(1R,2R)-2-hydroxy-1-(hydroxymethyl)propyl]-L-cysteinamide, cyclic (2→7)-disulfide, acetate

**MF:** C<sub>69</sub>H<sub>94</sub>N<sub>14</sub>O<sub>17</sub>S<sub>2</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

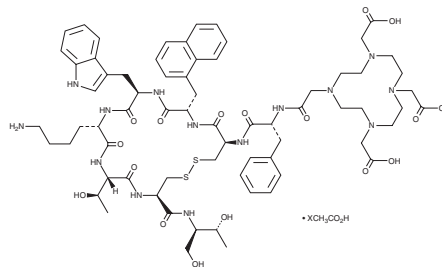
**FW:** 1,455.7

**Purity:** ≥98%

**Supplied as:** A 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

DOTA-NOC (acetate) is supplied as a solid. Aqueous solutions of DOTA-NOC (acetate) can be prepared by directly dissolving the solid in aqueous buffers. DOTA-NOC (acetate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

DOTA-NOC is a synthetic peptide composed of a derivative of the somatostatin receptor (SST) agonist octreotide (Item No. 23757) conjugated to the bifunctional chelator DOTA.<sup>1</sup> It has been conjugated to radioligands for use in PET imaging of neuroendocrine tumors and for tissue distribution studies *in vivo*.<sup>1,2</sup>

### References

1. Wild, D., Schmitt, J.S., Ginj, M., *et al.* DOTA-NOC, a high-affinity ligand of somatostatin receptor subtypes 2, 3 and 5 for labelling with various radiometals. *Eur. J. Nucl. Med. Mol. Imaging* **30(10)**, 1338-1347 (2003).
2. Dam, J.H., Langkjær, N., Baun, C., *et al.* Preparation and evaluation of [<sup>18</sup>F]AlF-NOTA-NOC for PET imaging of neuroendocrine tumors: Comparison to [<sup>68</sup>Ga]Ga-DOTA/NOTA-NOC. *Molecules* **27(20)**, 6818 (2022).

#### 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, 11/14/2024

#### 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