

PRODUCT INFORMATION

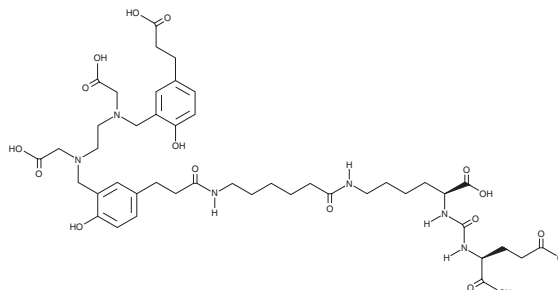


PSMA11

Item No. 42984

CAS Registry No.: 1366302-52-4
Formal Name: 22-[3-[[[2-[[[5-(2-carboxyethyl)-2-hydroxyphenyl]methyl](carboxymethyl)amino]ethyl](carboxymethyl)amino]methyl]-4-hydroxyphenyl]-5,13,20-trioxo-4,6,12,19-tetraazadocosane-1,3S,7S-tricarboxylic acid
Synonyms: Prostate-specific Membrane Antigen 11, PSMA-HBED-CC

MF: C₄₄H₆₂N₆O₁₇
FW: 947.0
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

PSMA11 is supplied as a solid. A stock solution may be made by dissolving the PSMA11 in the solvent of choice, which should be purged with an inert gas. PSMA11 is soluble (≥10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of PSMA11 can be prepared by directly dissolving the solid in aqueous buffers. PSMA11 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

PSMA11 is an unlabeled radiotracer specific for prostate-specific membrane antigen (PSMA) that is composed of a gallium chelator and a PSMA binding motif.¹ A radiolabeled form of PSMA11, [⁶⁸Ga]PSMA11, has been used as a PET tracer for *in vivo* imaging of PSMA⁺ tumors.

Reference

1. Eder, M., Schäfer, M., Bauder-Wüst, U., *et al.* ⁶⁸Ga-complex lipophilicity and the targeting property of a urea-based PSMA inhibitor for PET imaging. *Bioconjug. Chem.* **23**(4), 688-697 (2012)

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, 04/14/2025

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