

PRODUCT INFORMATION



Poly Citrulline (acetate)

Item No. 44481

Formal Name: (6S,9S,12S,15S,18S,21S,24S,27S,30S,33S)-1,6-diamino-1,7,10,13,16,19,22,25,28,31-decaoxo-9,12,15,18,21,24,27,30,33-nonakis(3-ureidopropyl)-2,8,11,14,17,20,23,26,29,32-decaazatetatriacontan-34-oic acid, acetate

Synonyms: Cit-Cit-Cit-Cit-Cit-Cit-Cit-Cit-OH, Decacitrulline

MF: $C_{60}H_{112}N_{30}O_{21} \cdot XC_2H_4O_2$

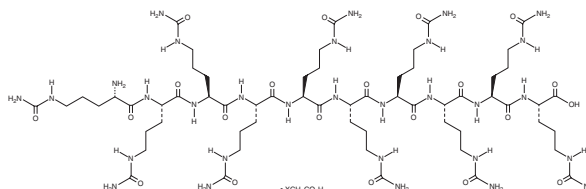
FW: 1,589.7

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Poly citrulline (acetate) is supplied as a solid. A stock solution may be made by dissolving the poly citrulline (acetate) in the solvent of choice, which should be purged with an inert gas. Poly citrulline (acetate) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Poly citrulline is a citrulline-containing decapeptide. It has been used as a hapten in the generation of anti-citrulline antibodies, which have been used to detect citrullination in postmortem human brain samples and synovial biopsy samples from patients with rheumatoid arthritis.^{1,2}

References

1. Makrygiannakis, D., Revu, S., Engström, M., *et al.* Local administration of glucocorticoids decreases synovial citrullination in rheumatoid arthritis. *Arthritis Res. Ther.* **14**(1), R20 (2012).
2. Nicholas, A.P. and Whitaker, J.N. Preparation of a monoclonal antibody to citrullinated epitopes: Its characterization and some applications to immunohistochemistry in human brain. *Glia* **37**(4), 328-336 (2002).

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.

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