

Ubiquilabel™, Ubiquitin [¹⁵N (98%), ²H (99%)]

Cat. # NS103

Background: ¹⁵N and ²H, heavy-labeled ubiquitin provides a stable standard for ubiquitin structural study by NMR. The recombinant human ubiquitin is purified from *E.coli* expressing the human ubiquitin gene. The cells are grown in media where the sole nitrogen source is ¹⁵N labeled ammonium sulfate and respective amount of ²H. The recombinant human ubiquitin is purified without heating by a proprietary procedure that preserves the native structure of the protein. The purified protein is homogenous when analyzed by overloaded SDS gels. The pools of purified protein solution are extensively dialyzed against water and lyophilized.

Application: Heavy-labeled ubiquitin for NMR studies

Product Information

Purity:	98% by RP-HPLC
Protein:	Ubiquitin
Atom% Enrichment:	¹⁵ N (98%), ² H (99%)
Physical State:	Lyophilized powder
Quantity:	5mg or 10mg
Solubility:	>30mg/ml in aqueous solution
Storage:	-20°C

References

1. Tjandra, N., Feller, S.E., Pastor, R.W. & Bax, A. Rotational diffusion anisotropy of human ubiquitin from ¹⁵N NMR relaxation. *J. Am. Chem. Soc.* **117**, 12562-12566 (1995)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved