

**Ubiquilabel™, Ubiquitin [<sup>15</sup>N (98%)]**

Cat. # NS101

**Background:** <sup>15</sup>N, 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 <sup>15</sup>N labeled ammonium sulfate. 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**

<b>Purity:</b>	98% by RP-HPLC
<b>Protein:</b>	Ubiquitin
<b>Atom% Enrichment:</b>	<sup>15</sup> N (98%)
<b>Molecular Weight:</b>	8,668.25 Da
<b>Physical State:</b>	Lyophilized powder
<b>Quantity:</b>	5mg or 10mg
<b>Solubility:</b>	>30mg/ml in aqueous solution
<b>Storage:</b>	-20°C

**References**

1. Tjandra, N., Feller, S.E., Pastor, R.W. & Bax, A. Rotational diffusion anisotropy of human ubiquitin from <sup>15</sup>N 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