

Single Lysine Ubiquitin Explorer Panel

Cat. # SI210

Background: Ubiquitin is a small polypeptide that can be conjugated via its C-terminus to the primary amine groups of a lysine residue on target proteins. This covalent attachment is referred to as monoubiquitylation. Additional ubiquitin moieties can subsequently be conjugated to this initial ubiquitin, utilizing any one (of seven) lysine residue on the surface of ubiquitin. The formation of these ubiquitin chains is referred to as polyubiquitylation.

LifeSensors has produced a complete panel of tag-free recombinant human ubiquitins that are engineered to have all available lysines mutated to arginines except at position mentioned. For example, ubiquitin-K6 has all available lysines mutated to arginines except at position 6, such that this molecule can only form polyubiquitin chains of K6-linkage type. The panel consists of 1 mg of each Ub mutant- K6, K11, K27, K29, K33, K48, and K63. The panel also contains 1 mg each of wt Ub and ubiquitin-K0 (all lysines mutated to arginines).

General Product Information

| | |
|--------------------|---|
| Purity: | ≥ 95% by SDS-PAGE |
| Quantity: | 1 mg each |
| Solubility: | >50mg/mL |
| Storage: | -80° C. Avoid repeated freeze/thaw cycles |

| Kit Contents | Amount | Buffer |
|---------------|--------|------------|
| Wt Ub | 1 mg | PBS pH 7.4 |
| Ubiquitin-K0 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K6 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K11 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K27 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K29 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K33 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K48 | 1 mg | PBS pH 7.4 |
| Ubiquitin-K63 | 1 mg | PBS pH 7.4 |

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2014 LifeSensors, Inc. All Rights Reserved