

# HEK293 cell lysate

Cell Lysate



Cat. No. 66-3002-010  
Lot. No. 30105

Quantity: 10 mg  
Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS Page 1 of 1

## Description

HEK293 cell lysate may be used to demonstrate the ability of ubiquitin binding domain proteins (Optineurin and NEMO) to capture polyubiquitylated proteins including IL-1 receptor-associated kinase 1 (IRAK1) (Windheim *et al.* 2008).

### Reference:

Windheim M, Stafford M, Peggie M, Cohen P (2008) Interleukin-1 (IL-1) induces the Lys63-linked polyubiquitination of IL-1 receptor-associated kinase 1 to facilitate NEMO binding and the activation of IκappaBalpha kinase. *Mol Cell Biol* 28 1783-91.

## Physical Characteristics

**Species:** human

**Source:** unstimulated embryonic kidney (HEK293) cells

**Quantity:** 10 mg

**Concentration:** 20 mg/ml

**Formulation:** 50 mM Tris/HCl pH7.5, 1 mM EGTA, 1 mM EDTA, 1% Triton X-100, 1 mM sodium orthovanadate, 50 mM sodium fluoride, 5mM sodium pyrophosphate, 10 mM sodium β-glycerophosphate, 270 mM sucrose, 1 mM PMSF, 1 mM benzamidine, 50 mM iodoacetamide

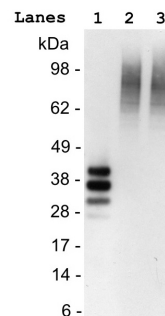
**Stability/Storage:** 12 months at -70°C; aliquot as required

## Quality Assurance

### Identification of Polyubiquitin Chains in Lysate:

A Western blot of IL-1 stimulated HEK293 cell lysate (lane 2) and non-stimulated HEK293 cell lysate (lane 3) probed with an anti-FK2 antibody indicates that both of the lysates are abundant with ubiquitylated proteins and possibly free ubiquitin chains. Polyubiquitin (Ub2-7) (K63-linked) was included as a positive control in this experiment (lane 1).

|                                     |   |   |   |
|-------------------------------------|---|---|---|
| Poly-ubiquitin (Ub2-7) (K63-linked) | + | - | - |
| HEK293 lysate (IL-1 stimulated)     | - | + | - |
| HEK293 lysate                       | - | - | + |



www.ubiquigent.com  
Dundee, Scotland, UK

### ORDERS / SALES SUPPORT

International: +1-617-245-0003  
US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233)  
Email: sales.support@ubiquigent.com

### UK HQ and TECHNICAL SUPPORT

International: +44 (0) 1382 381147 (9AM-5PM UTC)  
US/Canada: +1-617-245-0020 (9AM-5PM UTC)  
Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2012. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0