

# ULK1 Antibody FITC

## Ordering Information

[Order Online »](#)

|               |       |
|---------------|-------|
| FGX-ULK1-FITC | 100µg |
|---------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

|                           |   |
|---------------------------|---|
| <b>Application</b>        | ELISA, IF, IP, WB   |
| <b>Host</b>               | Rabbit  |
| <b>Immunogen</b>          | Synthetic peptide taken within amino acid region 1000-1050 on mouse ULK1 protein. |
| <b>Isotype</b>            | IgG   |
| <b>Purity Detail</b>      | Affinity Purified   |
| <b>Species Reactivity</b> | Bovine, Human, Monkey, Mouse, Panda   |
| <b>UniProt ID</b>         | O75385  |



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)