

Mono- and polyubiquitinated conjugates recombinant monoclonal antibody (UBC-JNK1/2/3/4/6/7/8/9/10/11/12/13/14/15/16/17/18/19/20/21/22/23/24/25/26/27/28/29/30/31/32/33/34/35/36/37/38/39/40/41/42/43/44/45/46/47/48/49/50/51/52/53/54/55/56/57/58/59/60/61/62/63/64/65/66/67/68/69/70/71/72/73/74/75/76/77/78/79/80/81/82/83/84/85/86/87/88/89/90/91/92/93/94/95/96/97/98/99/100) (UBC-JNK1/2/3/4/6/7/8/9/10/11/12/13/14/15/16/17/18/19/20/21/22/23/24/25/26/27/28/29/30/31/32/33/34/35/36/37/38/39/40/41/42/43/44/45/46/47/48/49/50/51/52/53/54/55/56/57/58/59/60/61/62/63/64/65/66/67/68/69/70/71/72/73/74/75/76/77/78/79/80/81/82/83/84/85/86/87/88/89/90/91/92/93/94/95/96/97/98/99/100) conjugate)

This antibody is covered by our [Warranty-Free Guarantee](#).

Citations: [1](#)

[View Online »](#)

[Order Online »](#)

ENZ-ABS840F-0025	25µg
ENZ-ABS840F-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Flow Cytometry, ICC
Clone	UBCJ2
Formulation	Liquid. In PBS containing 0.01% sodium azide.
Host	Mouse
Immunogen	Poly-ubiquitinated lysozyme.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:250)Flow cytometry (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from tissue culture supernatant.
Species Reactivity	Species independent
Specificity	Recognizes mono- and polyubiquitinated protein conjugates in a wide range of species.

UniProt ID P0CG47 (UBB), P0CG48 (UBC)



Worry-free Guarantee

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +32 3 466 0420
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

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

This antibody is covered by our Worry-Free Guarantee