



Mouse (monoclonal) anti-ERK5/MAPK7

PRODUCT ANALYSIS SHEET

Catalog Number:	44688M
Volume:	50 µL
Clone Number:	12F2
Isotype:	IgG1 (mouse)
Form of Antibody:	Mouse monoclonal immunoglobulin in PBS, pH 7.3, with PEG and Sucrose.
Preservative:	0.09% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Purification:	The antibody was purified from serum-free cell culture supernatant by subsequent ultrafiltration and size exclusion chromatography.
Immunogen:	Peptide (N-terminal sequence) conjugated to hemocyanin.
Target Summary:	Extracellular signal-regulated kinase 5 (ERK5, MAPK7, big mitogen activated kinase 1, BMK1) belongs to the serine/threonine protein kinase family. It contains a TEY motif in the activation loop similar to ERK1/2. Extracellular signals, including receptor tyrosine kinases and G-protein-coupled receptors as well as osmotic and oxidative stress lead to ERK5 activation by MEK5. The ERK5 pathway plays an important role in cellular proliferation, differentiation and survival
Specificity:	This Mab recognizes ERK5/MAPK7 (92 kDa).
Species Reactivity:	Human, mouse, rat and dog.
Applications:	The antibody is suitable for Western blotting. Other applications may be possible but have not been tested.
Suggested Working Dilutions:	Immunoblotting: 0.5 µg/mL for HRP/ECL detection. The optimal antibody concentration should be determined empirically for each specific application.
Recommended Positive Control:	Cell lysate from untreated SKOV-3 cells.
Storage:	Upon arrival, we recommend a brief centrifugation before opening to settle vial contents. Then, apportion the antibody into working aliquots and store at -20°C. Avoid repeated freeze / thaw cycles.
Expiration Date:	Expires one year from date of receipt when stored as instructed.

This product is for research use only. Not for use in diagnostic procedures.

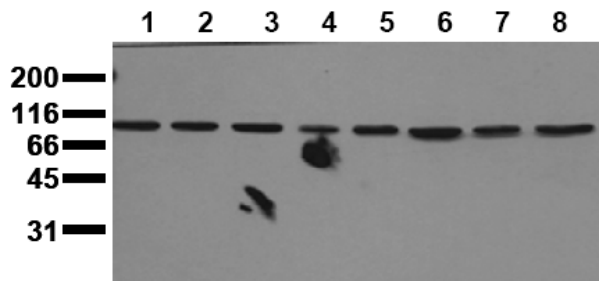
www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI44688M

(Rev 11/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



Western Blotting

Extracts of serum starved A431 (1), A549 (2), SKOV-3 (3), OVCAR-5 (4), HaCaT (5), PC3 (6), HeLa (7) and HepG2 (8) tumor cells (approximately 20,000 cells per lane) were resolved by SDS-PAGE and transferred to PVDF. The membrane was blocked with a casein/Tween 20 buffer then incubated with the Mab at 0.5 $\mu\text{g}/\text{mL}$ for 1 hour at room temperature. After washing, the membrane was incubated with an anti-mouse HRP-conjugated secondary antibody and signals were detected using an ECL detection method (exposure time: 30 seconds).

The data show that the Mab recognizes various levels of ERK5/MAPK7 in these cell systems.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI44688M

(Rev 11/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.