



Qty: 100 µg/400 µL

Rabbit anti-Connexin 31.1 (Mid)

Catalog No. 487800

Lot No.

## Rabbit anti-Connexin 31.1 (Mid)

### FORM

This polyclonal antibody is supplied as a 400 µL aliquot at a concentration of 0.25 mg/mL in phosphate buffered saline (pH 7.4) containing 0.1% sodium azide. This antibody is epitope-affinity purified from rabbit antiserum.

PAD: ZMD.693

### IMMUNOGEN

Synthetic peptide derived from the internal region of the mouse Connexin 31.1 protein (Accession# NP\_034421), which is 82% homologous to rat sequence.

### SPECIFICITY

This antibody is specific for the Connexin 31.1 (gap junction protein beta 5, Gjb5, Cx31.1, Cnx31.1) protein. On Western blots, it identifies the target band at ~31 kDa. The specificity of this antibody has been validated by tissue samples from the Connexin 31.1 knockout mice.

### REACTIVITY

Reactivity has been confirmed with Connexin 31.1 transfected HeLa cell lysates and mouse skin homogenates by Western blotting and with mouse skin frozen sections by immunohistochemistry. Skin tissue from the Connexin 31.1 knockout mice has been used as the negative control. Reactivity with rat is expected.

Sample	Western Blotting	Immuno-cytochemistry	Immuno-histochemistry
Human	ND	ND	ND
Mouse	+++	+++	++
Rat	ND	ND	ND

(Excellent +++, Good++, Poor +, No reactivity 0, Not applicable N/A, Not Determined ND)

### USAGE

Working concentrations for specific applications should be determined by the investigator. Appropriate concentrations will be affected by several factors, including secondary antibody affinity, antigen concentration, sensitivity of detection method, temperature and length of incubations, etc. The suitability of this antibody for applications other than those listed below has not been determined. The following concentration ranges are recommended starting points for this product.

**Western Blotting:** 2-4 µg/mL  
**Immunocytochemistry:** 2-4 µg/mL  
**Immunohistochemistry:** 2-4 µg/mL

### STORAGE

Store at 2-8°C for up to one month. Store at -20°C for long-term storage. Avoid repeated freezing and thawing.

(cont'd)

[www.invitrogen.com](http://www.invitrogen.com)

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

PI487800

(Rev 10/08) 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](http://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.

**BACKGROUND**

Gap junctions are conduits that allow the direct cell-to-cell passage of small cytoplasmic molecules, including ions, metabolic intermediates, and second messengers, and thereby mediate intercellular metabolic and electrical communication.<sup>1</sup> Gap junction channels consist of connexin protein subunits, which are encoded by a multigene family. Connexin 31.1 is a gap junction protein of 271 amino acids encoded by a novel single copy mouse gene on mouse chromosome 4.<sup>2</sup> Northern blot analysis revealed predominant expression of Connexin 31.1 in skin and two related mouse keratinocyte cells. Minor levels of Connexin 31.1 mRNA were detected in testis.<sup>2</sup> Connexin 31.1 mRNA was 15-30 times more abundant in mature skin than in fetal skin.<sup>3</sup> In response to epidermal wounding, Connexin 31.1 was down-regulated in cells both peripheral to and at the wounded edge. The data thus suggest that intercellular communication is intimately involved in regulating epidermal wound repair.<sup>4</sup>

**REFERENCES**

1. Nagy JI, et al. *Brain Res Brain Res Rev* 47(1-3):191-215, 2004.
2. Hennemann H, et al. *J Biol Chem* 267(24):17225-17233, 1992.
3. Goliger JA, et al. *Dev Dyn* 200(1):1-13, 1994.
4. Goliger JA, et al. *Mol Biol Cell* 6(11):1491-1501, 1995.

**RELATED PRODUCTS**

<b>Product</b>	<b>Conjugate</b>	<b>Cat. No.</b>
Protein A	Sepharose 4B	10-1041
rec-Protein G	Sepharose 4B	10-1241
ZyMAX™ Goat anti-rabbit IgG	Unconjugated	81-6100
ZyMAX™ Goat anti-mouse IgG	Unconjugated	81-6500

Secondary antibody conjugates.

<b>Conjugate</b>	<b>Goat anti-rabbit IgG (H+L)</b>	<b>Goat anti-mouse IgG (H+L)</b>	<b>Ex/Em*</b>	<b>Fluorescence similar to--</b>
Alexa Fluor® 488	A11008	A11001	495/519	FITC
Alexa Fluor® 555	A21428	A21422	555/565	Cy3
Alexa Fluor® 594	A11012	A11005	590/617	Texas Red
Alexa Fluor® 647	A21244	A21235	650/668	Cy5
HRP	81-6120	81-6520	NA**	NA
AP	81-6122	81-6522	NA	NA
Biotin	B2770	B2763	NA	NA

\*Excitation/emission (nm); \*\*Not applicable

For additional secondary antibody conjugates, visit [www.invitrogen.com/antibodies](http://www.invitrogen.com/antibodies)

**For Research Use Only**

[www.invitrogen.com](http://www.invitrogen.com)

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

PI487800

(Rev 10/08) 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](http://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.