

Fn14 (mouse):Fc (human), (recombinant)

The fibroblast growth factor-inducible-14 (Fn14) is a 129 aa type Ia transmembrane protein. Fn14 is a growth-factor regulated, immediate-early response protein. It is expressed in a developmental and adult tissue-specific manner. Fn14 expression decreases cellular adhesion to extracellular matrix proteins fibronectin and vitronectin and reduces cell growth and migration. It is expressed at high levels in liver cancer cells, in hepatocellular carcinoma specimens as well as in regenerating liver cells showing Fn14 may play a role in hepatocyte growth and liver neoplasia. Fn14 has been described to bind TWEAK (Apo-3 Ligand), a member of the TNF family. Historical data has shown that Fn14 inhibits TWEAK-mediated killing of Kym-1 target cells.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-036-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TWEAK-R, TNFRSF 12A, Fibroblast growth factor-inducible immediate-early response protein 14, CD266
Application Notes	ELISA: binds to TWEAK.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; BioWhittaker).
Formulation	Lyophilized. Contains PBS.
Gene/Protein Identifier	A003364 (UCSD Signaling Gateway ID)
MW	~34kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in HEK 293 cells. The extracellular domain of mouse Fn14 (aa 1-75) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse TWEAK.
UniProt ID	Q9CR75



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)