## IGF-I (human), (recombinant)

Recombinant human IGF-I that promotes cell proliferation and inhibits apoptosis.

Insulin-like growth factor-1 (IGF-1) was originally known as Somatomedin C but was later changed to IGF-1 due to its structural similarity to proinsulin. IGF-1 has insulinlike effects, but it appears to be involved in many other functions and has been linked to aging, nutrition, physical activity and cancer. While each of these is very interesting on its own, together they may shed light on the role that lifestyle plays in influencing cancer risk. High levels of circulating IGF-1 and low levels of IGF binding protein-3 (IGFBP-3) are associated with increased risk of several cancers. IGF-1 can stimulate cell proliferation and inhibit apoptosis in many differe nt kinds of cells. IGF-1 circulates in blood bound with IGFBPs. Six IGFBPs have been identified but the majority of circulating IGF -1 is complexed with IGFBP-3. The IGF-1 and IGFBP-3 exist in a complex with a third protein, named the acid -labile subunit.

Citations: 1

View Online »

**Ordering Information** 

**Order Online** »

ADI-908-059-0100

100µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** Lyophilized IGF-I although stable at room temperature for 3 weeks, should be stored

desiccated below -18°C. Upon reconstitution IGF-I should be stored at +4°C between 2-7 days and for future use below -18°C. For long term storage it is recommended to add a

carrier protein (0.1% HSA or BSA).

**Handling** Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Insulin-like growth factor I

Application Notes Western blot control.

Biological Activity

The activity as determined by the dose-dependent

proliferation of mouse FDC-P1 is less than 1.0 ng/ml, corresponding to a specific activity of 1×10<sup>6</sup> units/mg.

Formulation Lyophilized from a sterile filtered concentrated solution in

PBS, pH 7.0.

MW ~8kDa

Purity ≥95% (SDS and HPLC)

Purity Detail Purified by SDS and HPLC.

**Reconstitution** Reconstitute the lyophilized IGF-1 in sterile 18MΩ-cm H2O

not less than 100μg/ml, which can then be further diluted

to other aqueous solutions.

Source Produced in E. coli.



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com P05019
Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

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

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

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