

Human entry ORF set—cytokine

Cat. #OHS4929

Vector: pDONR223
Number of clones: 166
Number of genes: 130
Species: *Homo sapiens*
Ontology: Molecular function
GO term: Cytokine activity
GO ID: GO:0005125
Source: CGAP Gene Finder

Product description

This clone set contains Dharmacon Human Entry ORFs whose annotations fall within the Gene Ontology (GO) term indicated above. The source collection is the Dharmacon Human ORF Collection (Releases 1+2) created by the Dana-Farber Cancer Institute. In some cases, there will be multiple ORF clones for the same gene within the set. These usually correspond to alternatively spliced mRNA isoforms, but it is possible that a few are redundant. In other cases, ORFs may be derived from mRNA isoforms that lack the functional domain annotated for the corresponding gene locus. No effort has been made to remove redundant clones or clones lacking a specific functional domain. A complete clone list is contained in the accompanying data file.*

If you have any questions, contact

t +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f +44 (0)1223 655 581

w horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

About GO

GO provides structured controlled vocabularies (ontologies) to describe gene attributes in any organism and is a useful framework for assembly of cDNA clones into sets. The three ontologies (biological process, cellular component and molecular function) are each a hierarchy of terms organized from general to specific classifications. GO annotations are applied to genes by curators within the Gene Ontology Consortium and do not originate from Dharmacon, Inc. Consult the Gene Ontology website for details (geneontology.org).

Quality control

All cultures are checked for growth prior to shipment.

Shipping and storage

The set is shipped as bacterial cultures of *E. coli* glycerol stocks arrayed into 96-well plates. The set is shipped on dry ice and should be immediately stored at -80°C upon receipt.

To allow any CO_2 that may have dissolved into the media from the dry ice in shipping to dissipate, please store constructs at -80°C for at least 48 hours before thawing.

*Data file is available on the USB provided with product shipment and is also available at dharmacon.horizondiscovery.com.