

Human CD40L(CD154)-transfected murine L cells

Product reference: DDX-S2

huCD40L/huCD154 transfected murine L cells (L6)

Murine L cells were stably transfected with human CD40L (L6) and freeze-dried through proprietary process. These cells are primarily intended to activate human cells through triggering of CD40 surface antigen but were also demonstrated to activate B cells from several animal species. L6 cells successfully support the differentiation of human monocyte towards the dendritic cell lineage.

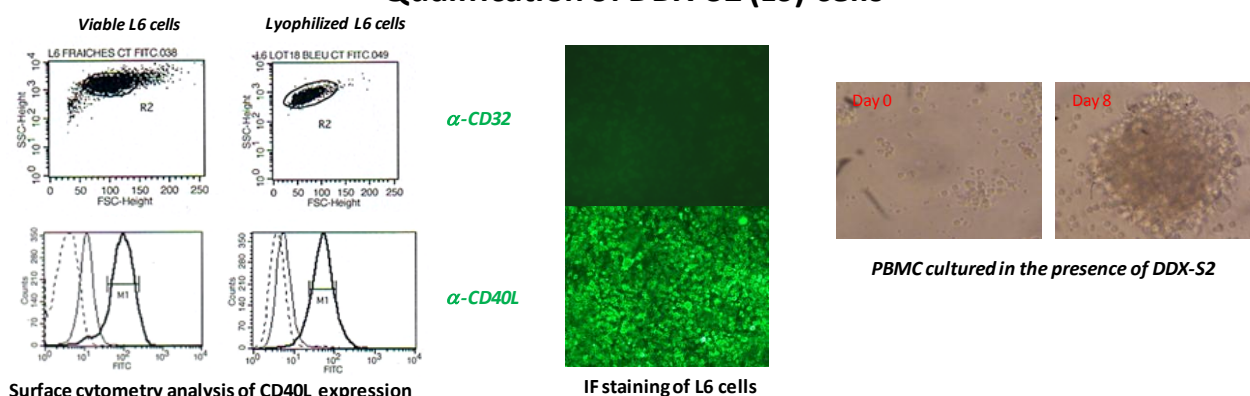
L6 cells are included in the Human Blood B booster and the Animal Blood B Booster kits, and support B cell growth and Ig secretion, but can also used alone for transient cultures.

Available formats:

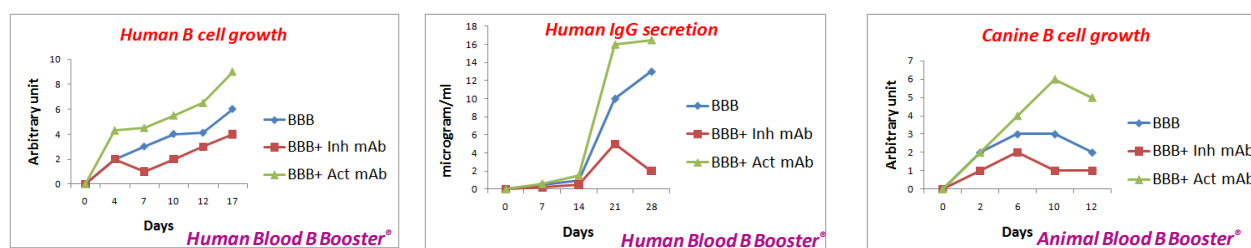
| Reference | Format | Size | Application |
|-----------|--------------|--------------------------|--|
| DDX-S2-2 | freeze-dried | 2.10 ⁶ cells | Activation of CD40 ⁺ human cells, in vitro generation of monocyte-derived dendritic |
| DDX-S2-5 | | 5.10 ⁶ cells | |
| DDX-S2-10 | | 10.10 ⁶ cells | |

Resuspended L6 cells can be kept at 4°C under sterile conditions for up to 2 months.

-Qualification of DDX-S2 (L6) cells

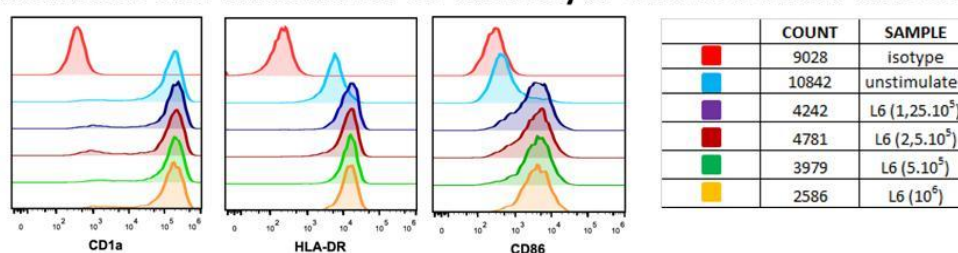


- DDX-S2 supports B cell growth and Ig secretion



Freeze-dried L6 cells are part of the Blood B Booster technology intended to immortalize human B cells and to activate animal B cells

-DDX-S2 promotes the maturation of monocyte-derived dendritic cells in vitro



Monocytes isolated from PBMC were cultured 5 days in the presence of rhuGM-CSF (200ng/ml) and rhuIL4 (10ng/ml). 10⁵ sorted CD1a⁺ cells were cultured for 24h with increasing number of L6 cells. Facs staining of CD1a, HLA-DR and CD86 are shown.

Storage conditions

-20°C. KEEP CONTENTS STERILE: no preservative.

Usage recommendation:

Resuspension in sterile complete culture medium

Not for use in Humans. For research purpose only