

Product	AbID	Standard ID	NIBSC conversion factor
Human Erythropoietin (EPO)	ab274397	11/170	0.001
Human AFP	ab193765	AFP	0.927
Human alpha 1 Antitrypsin (SERPINA1)	ab189579	05/162	0.3649
Human BMP2	ab277085	93/574	0.00024
Human C-Peptide	ab260064	13/146	1.077
Human Flt-3 ligand	ab254507	96/532	0.0014
Human GM-CSF	ab174448	88/646	0.0078
Human Growth Hormone	ab190811	80/505	2.3x10-6
Human HGF (Hepatocyte Growth Factor)	ab275901	96/564	0.001121
Human HIV1 p24	ab218268	90/636	0.43
Human IFN gamma High Sensitivity	ab236895	87/586	0.01113
Human IGFBP3	ab211652	93/560	0.828
Human IgM	ab214568	67/086	0.000249
Human IL-12 p70	ab223592	95/544	0.012
Human IL-15	ab218266	95/554	0.036
Human IL-17A	ab216167	01/420	0.0496
Human IL-2	ab270883	86/500	1.4
Human IL-29	ab236715	10/176	0.03255
Human IL-4	ab215089	88/656	0.0167
Human IL-5	ab215536	90/586	0.00613
Human IL-7	ab270218	90/530	0.266
Human IL-8	ab214030	89/250	0.0042
Human Interferon alpha 1	ab213479	94/784	(no conversion factor)
Human Interferon beta	ab278127	00/572	0.5342
Human Interferon omega	ab282309	94/754	0.1203
Human MCP1	ab179886	92/794	0.000266
Human M-CSF	ab245714	89/512	0.89
Human Thrombopoietin	ab219632	03/124	0.0322
Human TNF alpha	ab181421	12/154	0.094
Human Transferrin Receptor (CD71)	ab272469	07/202	0.01144
Human VEGF	ab222510	02/286	0.0011
Human Von Willebrand Factor	ab223864	07/316	0.3238
Mouse IL-2	ab223588	93/566	0.3051