

Currency conversions calculated December 30, 2021

By the numbers	Simple Plex Assay	Traditional ELISA
Cost per Assay	Euro (EUR): €698 Japanese yen (JPY): ¥90,958 British Pound sterling (GBP): £585	Euro (EUR): €485 Japanese yen (JPY): ¥63,332 British Pound sterling (GBP): £408
Format	72x1 cartridge	96-well plate
Wells for Standard Curve	0	16 (8-point curve in duplicate)
Available Wells for Samples	72	80
Number of Reportable Results	72 (each sample run in triplicate)	40 (each sample run in duplicate)
Cost per Reportable Result	Euro (EUR): €10 Japanese yen (JPY): ¥1,263 British Pound sterling (GBP): £8	Euro (EUR): €12 Japanese yen (JPY): ¥1,583 British Pound sterling (GBP): £10

Not seeing your currency conversion? Or want to learn more about the cost analysis?

[Request Simple Plex Pricing](#)