

HTRF[®] readout - Set up recommendations for Artemis readers

Two sequential measurements will be carried out: at 620 nm for the cryptate emission, and at 665 nm for the specific signal emitted by the acceptor (XL665 or d2). A ratio of the two fluorescence intensities* (acceptor/donor) then allows the calculation of Delta F (%), i.e. the relative energy transfer rate for each data point.

Artemis readers must be appropriately configured for HTRF[®] readout by setting up the measurement conditions in the "test protocol setting" function of the software piloting the reader. These parameters should be specifically entered as defined in the table below. No special upgrade is required for HTRF[®] readout, as it is a dedicated instrument.

Delay time	100 μ s
Integration time	100 μ s
Flash number	50

** The fluorescence ratio is a correction method developed by Cisbio international with an application limited to the use of HTRF[®] reagents and technology, and for which Cisbio international has granted a licence to Furuno. The method is covered by the US patent 5,527,684 and its foreign equivalents.*