

[E297] Extracellular OCR Plate Assay Kit

Please check the following notice before purchase and use!

1. About the Plate Reader

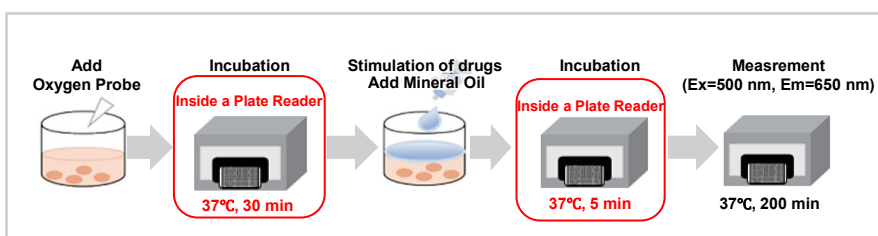
Please use a plate reader with the following functions:

- Temperature control
- Ex=500 nm, Em=650 nm filter
- Bottom Reading

2. Precautions

Please incubate inside a plate reader. If incubate with a incubator, the temperature difference may affect the OCR results and decrease the data reproducibility. For details of protocol, please refer to the product manual.

General Protocol



Please refer to the Step 3,7 with floating cells, Step 5,9 with adherent cell in the manual.

Incubate with a incubator instead of a plate reader will affect the result.

