

XF Glycolysis Stress Test with Astrocyte cells

ASSAY OVERVIEW: General guidelines for performing the **XF Glycolysis Stress Test** assay with Astrocyte cells. This **Assay Guide** is for use on either XF^e96, XF96, or XFp Analyzers[‡].

- This guide is associated with the **XF^e96** Assay Template: **Astrocyte-GLYCO-96** (.asyt file).
- This assay may be adapted for acute injections (compounds). Assign the acute injection to Port A, reassign the injections of glucose, oligomycin and 2-deoxyglucose to ports B, C and D, respectively.
- Cells are to be plated at the indicated density 1-2 day(s) prior to the assay; **primary cells are to be cultured for 10-14 days in separate vessel, then seeded in XF Cell Culture Microplate(s).*
- The compound concentrations listed are *final* concentrations in well.
- Sample data is provided below. Absolute rates and magnitude of responses may vary based on biological and experimental variables.

Please note: Further optimization may be required depending on parameters tested and variables modified.

INJECTION STRATEGY: XF Glycolysis Stress Test
(Final concentration in well)

- Port A: 10 mM glucose
- Port B: 1.0 μ M oligomycin
- Port C: 50 mM 2-deoxyglucose (2-DG)
- Port D: N/A

PRETREATMENTS:

- Control Group(s)
- Experimental Group(s)

ASSAY MEDIA: Glyco Stress Test Assay Medium

- XF Base Medium: Supplement with 2 mM glutamine, pH 7.4.
- Initial Assay Volume: 180-200 μ L

CELL SEEDING DENSITY:

- Astrocyte cells.
- 2.0×10^4 cells/well, plated 1-2 days prior to assay*.

INSTRUMENT PROTOCOL:

- Calibrate
- Equilibrate
- Basal: 3 cycles
- 3 min *Mix*, 0 min *Wait*, 3 min *Measure*
- Inject Port A followed by 3 cycles
- 3 min *Mix*, 0 min *Wait*, 3 min *Measure*
- Inject Port B followed by 3 cycles
- 3 min *Mix*, 0 min *Wait*, 3 min *Measure*
- Inject Port C followed by 3 cycles
- 3 min *Mix*, 0 min *Wait*, 3 min *Measure*

XFp ANALYZER:

- All assay parameters (assay volumes, cell seeding density and all concentrations of media components and XF Cell Mito Stress Test compounds) remain unchanged.
- Groups are limited to 2 per plate (3 wells per group).

TYPICAL ASSAY DATA RESULTS FOR XF^e96, XF96, and XFp Analyzers (Prior to normalization)

Expected range of initial rate 11-17 mPH/min	Glucose response of initial rate*	Oligomycin response of initial rate*	2-Deoxyglucose response of initial rate*
	175%	375%	10%

**The indicated values represent a percentage of the initial rate and may vary +/-20%*

[‡] For XF^e24 and XF24 Analyzers, refer to **Assay Tech Hints: Modifying XF^e96 Parameters for XF^e24 and XF24 Analyzers**