

## XF Glycolysis Stress Test with Neo-Jurkat Cells

**ASSAY OVERVIEW:** General guidelines for performing the **XF Glycolysis Stress Test** assay with Neo-Jurkat cells. This **Assay Guide** is for use on either XF<sup>e</sup>96, XF96, or XFp Analyzers<sup>†</sup>.

- This guide is associated with the **XF<sup>e</sup>96 Assay Template: Neo-Jurkat-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 via centrifugation just prior to the assay in XF Glyco Stress Medium.
- 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 μM 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 μL

**CELL SEEDING DENSITY:**

- Neo-Jurkat cells.
- 1.0 x 10<sup>5</sup> cells/well, plated just prior to assay via centrifugation (250 x g, 1-2 minutes).

**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<sup>e</sup>96, XF96, and XFp Analyzers** (Prior to normalization)

Expected range of initial rate	Glucose response of initial rate*	Oligomycin response of initial rate*	2-Deoxyglucose response of initial rate*
16-24 mpH/min	500%	620%	20%

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

<sup>†</sup> For XF<sup>e</sup>24 and XF24 Analyzers, refer to **Assay Tech Hints: Modifying XF<sup>e</sup>96 Parameters for XF<sup>e</sup>24 and XF24 Analyzers**