

# Scoring Sheet **JBScreen Pentaerythritol**



Name: \_\_\_\_\_ Sample Name: \_\_\_\_\_ Temperature: \_\_\_\_\_

Plate ID: \_\_\_\_\_ Sample Concentration: \_\_\_\_\_ Reservoir Volume: \_\_\_\_\_

Set-up Date: \_\_\_\_\_ Sample Buffer : \_\_\_\_\_ Total Drop Volume: \_\_\_\_\_  $\mu$ l

Observation Date: \_\_\_\_\_ Sample : \_\_\_\_\_  $\mu$ l Reservoir : \_\_\_\_\_  $\mu$ l Additive: \_\_\_\_\_  $\mu$ l

	100 mM NaOAc pH 4,6			100 mM MES pH 6,5			100 mM HEPES pH 7,5			100 mM Tris pH 8,5			
	1	2	3	4	5	6	7	8	9	10	11	12	
A	25% PEP 426 100 mM NaOAc	35% PEP 426 100 mM NaOAc	45% PEP 426 100 mM NaOAc	25% PEP 426 100 mM MES	35% PEP 426 100 mM MES	45% PEP 426 100 mM MES	25% PEP 426 100 mM HEPES	35% PEP 426 100 mM HEPES	45% PEP 426 100 mM HEPES	25% PEP 426 100 mM Tris	35% PEP 426 100 mM Tris	45% PEP 426 100 mM Tris	PEP 426
B	25% PEP 426 100 mM NaOAc 50 mM MgCl <sub>2</sub>	35% PEP 426 100 mM NaOAc 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 426 100 mM NaOAc 400 mM KCl	25% PEP 426 100 mM MES 50 mM MgCl <sub>2</sub>	35% PEP 426 100 mM MES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 426 100 mM MES 400 mM KCl	25% PEP 426 100 mM HEPES 50 mM MgCl <sub>2</sub>	35% PEP 426 100 mM HEPES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 426 100 mM HEPES 400 mM KCl	25% PEP 426 100 mM Tris 50 mM MgCl <sub>2</sub>	35% PEP 426 100 mM Tris 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 426 100 mM Tris 400 mM KCl	
C	25% PEP 629 100 mM NaOAc	35% PEP 629 100 mM NaOAc	45% PEP 629 100 mM NaOAc	25% PEP 629 100 mM MES	35% PEP 629 100 mM MES	45% PEP 629 100 mM MES	25% PEP 629 100 mM HEPES	35% PEP 629 100 mM HEPES	45% PEP 629 100 mM HEPES	25% PEP 629 100 mM Tris	35% PEP 629 100 mM Tris	45% PEP 629 100 mM Tris	PEP 629
D	25% PEP 629 100 mM NaOAc 50 mM MgCl <sub>2</sub>	35% PEP 629 100 mM NaOAc 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 629 100 mM NaOAc 300 mM KCl	25% PEP 629 100 mM MES 50 mM MgCl <sub>2</sub>	35% PEP 629 100 mM MES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 629 100 mM MES 300 mM KCl	25% PEP 629 100 mM HEPES 50 mM MgCl <sub>2</sub>	35% PEP 629 100 mM HEPES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 629 100 mM HEPES 300 mM KCl	25% PEP 629 100 mM Tris 50 mM MgCl <sub>2</sub>	35% PEP 629 100 mM Tris 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEP 629 100 mM Tris 300 mM KCl	
E	25% PEE 270 100 mM NaOAc	35% PEE 270 100 mM NaOAc	45% PEE 270 100 mM NaOAc	25% PEE 270 100 mM MES	35% PEE 270 100 mM MES	45% PEE 270 100 mM MES	25% PEE 270 100 mM HEPES	35% PEE 270 100 mM HEPES	45% PEE 270 100 mM HEPES	25% PEE 270 100 mM Tris	35% PEE 270 100 mM Tris	45% PEE 270 100 mM Tris	PEE 270
F	25% PEE 270 100 mM NaOA 50 mM MgCl <sub>2</sub>	35% PEE 270 100 mM NaOAc 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 270 100 mM NaOAc 400 mM KCl	25% PEE 270 100 mM MES 50 mM MgCl <sub>2</sub>	35% PEE 270 100 mM MES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 270 100 mM MES 400 mM KCl	25% PEE 270 100 mM HEPES 50 mM MgCl <sub>2</sub>	35% PEE 270 100 mM HEPES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 270 100 mM HEPES 400 mM KCl	25% PEE 270 100 mM Tris 50 mM MgCl <sub>2</sub>	35% PEE 270 100 mM Tris 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 270 100 mM Tris 400 mM KCl	
G	25% PEE 797 100 mM NaOAc	35% PEE 797 100 mM NaOAc	45% PEE 797 100 mM NaOAc	25% PEE 797 100 mM MES	35% PEE 797 100 mM MES	45% PEE 797 100 mM MES	25% PEE 797 100 mM HEPES	35% PEE 797 100 mM HEPES	45% PEE 797 100 mM HEPES	25% PEE 797 100 mM Tris	35% PEE 797 100 mM Tris	45% PEE 797 100 mM Tris	PEE 797
H	25% PEE 797 100 mM NaOAc 50 mM MgCl <sub>2</sub>	35% PEE 797 100 mM NaOAc 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 797 100 mM NaOAc 400 mM KCl	25% PEE 797 100 mM MES 50 mM MgCl <sub>2</sub>	35% PEE 797 100 mM MES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 797 100 mM MES 400 mM KCl	25% PEE 797 100 mM HEPES 50 mM MgCl <sub>2</sub>	35% PEE 797 100 mM HEPES 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 797 100 mM HEPES 400 mM KCl	25% PEE 797 100 mM Tris 50 mM MgCl <sub>2</sub>	35% PEE 797 100 mM Tris 200 mM (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	45% PEE 797 100 mM Tris 400 mM KCl	

For questions and inquires please contact [xtals@jenabioscience.com](mailto:xtals@jenabioscience.com)