

Plate Record Template

Name of experiment:

Date of experiment:

	H	G	F	E	D	C	B	A	
1									
2									
3									
4									
5									
6									
7									
8									
9									
10									
11									
12									

Technical Assistance: call either 303-322-2254 or e-mail tservice@cytoskeleton.com