

表 グッド緩衝剤の利用最適pH範囲

pKa(20°C)	化合物	5	6	7	8	9	10	11
6.15	MES		5.5	7.0				
6.46	Bis-Tris		5.7	7.3				
6.60	ADA		5.8	7.4				
6.80	PIPES		6.1	7.5				
6.90	ACES		6.0	7.5				
6.95	MOPSO		6.2	7.4				
7.15	BES		6.6	8.0				
7.20	MOPS		6.5	7.9				
7.50	TES		6.8	8.2				
7.55	HEPES		6.8	8.2				
7.7	TAPSO		7.0	8.2				
7.85	POPSO		7.2	8.5				
7.9	HEPPSO		7.4	8.6				
8.0	EPPS		7.5	8.5				
8.15	Tricine		7.8	8.8				
8.35	Bicine		7.7	9.1				
8.4	TAPS		7.7	9.1				
9.5	CHES					8.6	10.0	
10.40	CAPS						9.7	11.1