

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