



表1 酸解離定数

pKa	EDTA	NTA	CyDTA	DTPA	EDTA-OH	GEDTA	TTHA*	DHEG (Bicine)	IDA	EDDA	DPTA-OH	NTP	Methyl-EDTA	HIDA	EDDP	EDTPO	NTPPO	BAPTA
a ₁	1.99	1.89	2.43	2.14	2.51	2.00	2.42	8.14	2.58	6.59	1.60	3.62	2.60	2.25	6.87	1.46	0.30	—
a ₂	2.67	2.49	3.52	2.38	5.31	2.68	2.95	—	9.33	9.58	2.60	4.24	3.03	8.75	9.60	2.72	1.50	—
a ₃	6.16	9.73	6.12	4.26	9.86	8.85	4.16	—	—	6.96	6.96	9.30	6.20	—	—	5.05	4.64	5.47
a ₄	10.26	—	11.70	8.60	—	9.46	6.16	—	—	9.49	—	—	10.84	—	—	6.18	5.86	6.36
a ₅	—	—	—	10.53	—	—	9.40	—	—	—	—	—	—	—	—	6.63	7.30	—
a ₆	—	—	—	—	—	—	10.19	—	—	—	—	—	—	—	—	7.43	12.10	—
a ₇	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	9.22	—	—
a ₈	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	10.60	—	—

表2 キレート安定度定数 (生成定数) logK_{MHL}

金属	EDTA	NTA	CyDTA	DTPA	EDTA-OH	GEDTA	TTHA*	DHEG (Bicine)	IDA	EDDA	DPTA-OH	NTP	Methyl-EDTA	HIDA	EDDP	EDTPO	NTPPO	BAPTA	CN
Ag ⁺	7.32	5.16	8.15	8.70	6.71	6.88	8.67 (13.89)	—	8.16	—	—	—	—	—	—	—	—	—	18.92
Al ³⁺	16.13	9.5	18.63	18.4	12.43	13.90	19.7 (28.9)	—	6.93	—	—	—	—	—	—	—	—	—	—
Am ³⁺	18.16	—	18.79	22.92	—	—	—	—	1.67	—	4.91	—	8.1	9.75	—	22.47	—	—	—
Ba ²⁺	7.76	4.83	8.64	8.63	5.54	8.41	8.22 (11.63)	—	—	—	—	—	—	3.42	—	11.14	—	5.77	—
Be ²⁺	9.27	7.11	10.81	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Bi ³⁺	27.9	—	31.2	29.7	21.8	23.8	—	—	—	—	—	—	—	—	—	—	—	—	—
Bk ³⁺	—	—	19.60	22.79	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Ca ²⁺	10.96	6.41	12.50	10.74	8.14	11.00	10.06 (14.16)	—	2.59	—	6.69	—	10.40	5.3	~1	6.93	6.68	6.97	—
Cd ²⁺	16.46	9.54	19.23	19.31	13.6	16.70	18.65 (26.85)	6.3	5.73	10.8	12.10	3.4	16.00	7.41	5.6	16.53	—	12.2	16.85
Ce ³⁺	15.98	10.83	16.76	20.50	14.11	15.70	19.20	7.5	6.18	7.48	—	—	16.79	8.50	—	21.10	—	—	—
Ce ⁴⁺	24.20	10.97	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Cl ³⁺	19.09	—	19.42	22.57	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Co ³⁺	18.45	—	18.81	22.99	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Co ²⁺	16.31	10.38	18.92	18.4	14.4	12.50	17.1 (28.8)	6.1	6.95	11.25	13.92	4.8	14.4	9.27	7.3	21.89	—	8.70	19.90
Co ³⁺	40.60	—	—	—	—	—	—	—	29.6	—	—	—	—	9.0	—	15.49	—	—	64.0
Cr ²⁺	13.61	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Cr ³⁺	23.40	>10	—	—	—	2.54	—	—	—	—	—	—	—	—	—	—	—	—	—
Cs ⁺	0.15	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Cu ²⁺	18.80	12.96	21.30	21.53	17.35	17.8	19.2 (32.6)	8.6	10.63	16.2	17.21	9.1	17.7	13.38	15.10	18.95	—	11.7	25.0
Dy ³⁺	18.30	11.74	19.69	22.82	15.30	17.42	23.29	7.6	6.88	8.31	—	—	19.09	8.88	—	—	—	—	—
Er ³⁺	18.38	12.03	20.20	22.74	15.42	17.40	23.19 (26.92)	7.7	7.09	8.59	—	—	19.61	9.24	—	—	—	—	—
Eu ³⁺	17.35	11.52	18.77	22.39	15.35	17.10	23.85	8.0	6.73	8.38	—	—	18.26	8.99	—	—	—	—	—
Fe ²⁺	14.33	8.84	16.27	16.55	12.2	11.92	17.1	4.31	5.80	—	12.90	—	14.30	6.78	6.3	—	—	—	24.0
Fe ³⁺	25.1	15.87	28.05	28.6	19.8	20.5	26.8 (40.5)	—	10.42	—	—	—	—	11.64	13.1	19.6	—	—	31.0
Fm ³⁺	—	—	—	22.7	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Ga ³⁺	20.27	13.6	22.91	23.0	16.9	—	—	—	—	—	—	—	—	9.01	—	—	—	—	—
Gd ³⁺	17.0	11.54	18.80	22.46	15.22	16.94	23.83	7.70	6.68	8.13	—	—	18.21	14.63	—	21.80	—	—	—
Hf ⁴⁺	29.5	20.34	—	35.40	—	—	19.08	—	—	—	—	—	—	—	—	—	—	—	—
Hg ²⁺	21.8	14.6	24.30	27.0	20.1	23.12	26.8 (39.1)	14.17	11.76	—	18.4	—	—	5.48	—	—	—	—	41.4
Ho ³⁺	18.05	11.90	19.89	22.78	15.32	17.38	23.59	7.6	6.97	8.42	—	—	19.30	9.14	—	21.85	—	—	—
In ³⁺	24.95	16.9	28.74	29.0	17.16	—	—	—	9.54	—	—	—	—	11.00	—	—	—	—	—
La ³⁺	15.50	10.36	16.75	19.48	13.46	15.79	22.22 (25.62)	7.3	5.88	7.04	—	—	16.42	8.00	5.8	20.15	—	—	—
Li ⁺	2.79	2.51	4.13	—	—	1.17	—	—	—	—	—	—	—	—	—	—	—	—	—
Ld ³⁺	19.65	12.49	20.91	22.44	15.88	17.81	—	7.5	7.61	9.09	—	—	20.56	9.50	—	—	—	—	—

* ()内は1:2キレート



*()内は1:2キレート

金属	EDTA	NTA	CyDTA	DTPA	EDTA-OH	GEDTA	TTHA*	DHEG (Bicine)	IDA	EDDA	DPTA-OH	NTP	Methyl-EDTA	HIDA	EDDP	EDTPO	NTPO	BAPTA	CN
Mg ²⁺	8.69	5.46	10.32	9.3	7.0	5.21	8.10(14.38)	1.15	2.94	3.90	5.30	<1	8.8	3.5	1.6	5.69	6.49	1.77	-
Mn ²⁺	14.04	7.44	16.78	15.60	10.7	12.3	14.68(21.19)	3.9	-	7.05	9.06	-	14.5	6.4	3.4	12.70	-	8.73	-
Mn ³⁺	24.8	-	28.9	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Mo ⁵⁺	6.36	-	17.68	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Mo ⁶⁺	19.5	-	-	-	-	1.38	-	-	-	-	-	-	-	-	-	-	-	-	-
Na ⁺	1.66	2.15	2.70	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Nb ⁵⁺	40.87	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Nd ³⁺	16.61	11.26	17.69	21.60	14.86	16.28	22.82(26.75)	7.6	6.50	8.30	-	-	17.54	8.80	-	21.47	-	-	-
Ni ²⁺	18.62	11.54	19.40	20.32	17.00	13.60	18.1(32.4)	7.7	8.9	-	16.63	5.8	14.2	10.2	9.3	15.30	-	7.79	22
NpO ₂ ²⁺	9.70	-	-	-	-	-	-	-	6.27	-	-	-	6.08	-	-	-	-	-	-
Pa ⁵⁺	8.19~	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Pb ²⁺	11.96	11.39	19.68	18.80	15.50	14.71	17.1(28.1)	7.5	7.45	-	11.4	-	17.3	9.41	-	-	-	11.3	10.3
Pd ²⁺	18.04	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Pd ⁴⁺	26.40	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Pm ³⁺	16.75	-	18.71	-	-	-	-	-	-	-	-	-	-	8.97	-	-	-	-	-
Pr ³⁺	16.40	11.07	17.23	21.07	14.61	16.05	23.45	7.7	6.44	7.84	-	-	17.17	8.64	-	-	-	-	-
Pu ³⁺	18.12	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Pu ⁴⁺	17.66	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Pu ⁶⁺	16.39	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Sb ³⁺	24.8	-	-	-	20.2	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Sc ³⁺	23.1	12.7	25.4	-	-	25.4	-	8.0	-	-	-	-	-	-	-	-	-	-	-
Sm ³⁺	16.7	11.53	18.63	22.34	15.28	23.85	23.81	7.8	6.64	8.28	-	-	17.97	9.10	-	22.39	-	-	-
Sn ²⁺	18.3	-	-	-	-	23.85	-	-	-	-	15.2	-	15.1	-	-	-	-	-	-
Sr ²⁺	8.6	4.98	10.54	9.68	6.92	8.50	9.26(12.70)	-	2.23	-	5.33	-	10.7	3.77	-	5.41	-	-	-
Tb ³⁺	17.81	11.59	19.30	22.71	15.32	17.27	23.61	8.0	6.78	8.18	-	-	18.64	9.08	-	-	-	5.13	-
Th ⁴⁺	23.2	12.4	29.25	28.78	18.5	-	31.9	7.8	9.32	-	-	-	-	10.7	-	-	-	-	-
Ti ³⁺	17.3	-	-	-	-	-	-	-	-	8.75	-	-	-	-	-	-	-	-	-
TiO ₂ ²⁺	17.3	12.3	19.9	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Ti(H ₂ O) ₆ ³⁺	20.43	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Tl ⁺	6.53	4.75	5.33	5.97	5.45	-	-	-	1.32	-	4.4	-	-	-	-	-	-	-	-
Tl ³⁺	22.5	18.0	38.3	48.0	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Tm ³⁺	18.62	12.20	20.46	22.72	15.59	17.48	-	-	7.22	-	-	-	20.08	9.35	-	21.41	-	-	35
U ⁴⁺	25.8	-	26.9	-	-	-	-	-	8.93	11.41	-	-	-	-	-	-	-	-	-
UO ₂ ²⁺	10.4	9.56	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
V ²⁺	12.70	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
VO ₂ ²⁺	18.77	-	-	-	-	-	-	-	9.01	-	-	-	-	-	-	-	-	-	-
V ⁵⁺	18.05	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
Y ³⁺	18.09	11.48	19.41	22.05	14.65	16.82	-	7.2	6.78	7.78	-	-	18.78	9.22	-	15.06	-	-	-
Yb ³⁺	18.88	12.40	20.80	22.62	15.88	17.78	23.58	7.7	7.42	8.93	-	-	20.25	9.38	-	-	-	-	-
Zn ²⁺	16.50	10.66	18.67	18.75	14.5	14.5	16.65(28.7)	6.5	7.27	11.22	13.70	5.3	16.2	8.33	7.6	17.05	-	9.38	16.9
Zr ⁴⁺	29.9	20.8	20.74	36.9	-	-	19.74	-	-	-	-	-	-	-	-	-	-	-	-

- キット類
- 酸化ストレス分子生物学
- 細胞内蛍光プローブ
- 細胞染色
- 菌染色用色素
- 細胞増殖/毒性
- 膜タンパク質
- ラベル化剤
- 二価性試薬
- 酸化還元
- イオン電極
- シンチレーター
- 生化学用緩衝剤
- キレート
- 比色/金属試薬
- 溶媒抽出
- 高純度溶媒
- その他
- SAM
- 機能的有機材料

*表示している希望納入価格は「本体価格のみ」で消費税等は含まれておりません。社会状況の変動により、予告なしに変更することがありますので、最新の価格はHPにてご確認ください。