



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	NTPO	BAPTA
a1	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	—
a2	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	—
a3	6.16	9.73	6.12	4.26	9.86	8.85	4.16	—	—	—	6.96	9.30	6.20	—	—	5.05	4.64	5.47
a4	10.26	—	11.70	8.60	—	9.46	6.16	—	—	—	9.49	—	10.84	—	—	6.18	5.86	6.36
a5	—	—	—	10.53	—	—	9.40	—	—	—	—	—	—	—	—	6.63	7.30	—
a6	—	—	—	—	—	—	10.19	—	—	—	—	—	—	—	—	7.43	12.10	—
a7	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	9.22	—	—
a8	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	10.60	—	—

表.2 キレート安定度定数 (生成定数) logK<sub>ML</sub>

\*( )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
Ag <sup>+</sup>	7.32	5.16	8.15	8.70	6.71	6.88	8.67(13.89)	—	—	—	—	—	—	—	—	—	—	—	18.92
Al <sup>3+</sup>	16.13	9.5	18.63	18.4	12.43	13.90	19.7(28.9)	—	8.16	—	14.4	—	—	—	—	—	—	—	—
Am <sup>3+</sup>	18.16	—	18.79	22.92	—	—	—	—	6.93	—	—	—	—	9.75	—	22.47	—	—	—
Ba <sup>2+</sup>	7.76	4.83	8.64	8.63	5.54	8.41	8.22(11.63)	—	1.67	—	4.91	—	8.1	3.42	—	11.14	—	5.77	—
Be <sup>2+</sup>	9.27	7.11	10.81	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Bi <sup>3+</sup>	27.9	—	31.2	29.7	21.8	23.8	—	—	—	—	—	—	—	—	—	—	—	—	—
Bk <sup>3+</sup>	—	—	19.60	22.79	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Ca <sup>2+</sup>	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 <sup>2+</sup>	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 <sup>3+</sup>	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 <sup>4+</sup>	24.20	10.97	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Cf <sup>3+</sup>	19.09	—	19.42	22.57	—	—	—	—	—	—	—	—	—	9.61	—	—	—	—	—
Cm <sup>3+</sup>	18.45	—	18.81	22.99	—	—	—	—	—	—	—	—	—	9.27	—	21.89	—	—	—
Co <sup>2+</sup>	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	9.0	7.3	15.49	—	8.70	19.90
Co <sup>3+</sup>	40.60	—	—	—	—	—	—	—	29.6	—	—	—	—	—	—	—	—	—	64.0
Cr <sup>2+</sup>	13.61	—	—	—	—	—	—	—	—	—	—	—	—	7.73	—	—	—	—	—
Cr <sup>3+</sup>	23.40	>10	—	—	—	2.54	—	—	—	—	—	—	—	—	—	—	—	—	—
Cs <sup>+</sup>	0.15	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Cu <sup>2+</sup>	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 <sup>3+</sup>	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 <sup>3+</sup>	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 <sup>3+</sup>	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 <sup>2+</sup>	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 <sup>3+</sup>	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 <sup>3+</sup>	—	—	—	22.7	—	—	—	—	—	—	—	—	—	—	—	—	—	—	—
Ga <sup>3+</sup>	20.27	13.6	22.91	23.0	16.9	—	—	—	—	—	—	—	—	9.01	—	—	—	—	—
Gd <sup>3+</sup>	17.0	11.54	18.80	22.46	15.22	16.94	23.83	7.70	6.68	8.13	—	—	18.21	—	—	21.80	—	—	—
Hf <sup>4+</sup>	29.5	20.34	—	35.40	—	—	19.08	—	—	—	—	—	—	14.63	—	—	—	—	—
Hg <sup>2+</sup>	21.8	14.6	24.30	27.0	20.1	23.12	26.8(39.1)	14.17	11.76	—	18.4	—	21.7	5.48	—	—	—	—	41.4
Ho <sup>3+</sup>	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 <sup>3+</sup>	24.95	16.9	28.74	29.0	17.16	—	—	—	9.54	—	—	—	—	11.00	—	—	—	—	—
La <sup>3+</sup>	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 <sup>+</sup>	2.79	2.51	4.13	—	—	1.17	—	—	—	—	—	—	—	—	—	—	—	—	—
Lu <sup>3+</sup>	19.65	12.49	20.91	22.44	15.88	17.81	—	7.5	7.61	9.09	—	—	20.56	9.50	—	—	—	—	—

