

Competitor Cross Guide:
Ellipsis vs. Cubis® II



Radwag Ellipsis

Features and Selling Points

- Limited experience in the weighing market
- Two year warranty
- 10" Touch Display
- Automatic sliding doors (IR Sensors from Semi-Micro upwards)
- Leveling:
 - Fully auto: from Semi-Micro upwards
 - (Assisted leveling) Semi-auto: Semi-Micro upwards
- Ionizer
- Internal two point adjustment
- Tool-free disassembly of chamber

User Log-In

- Password
- Camera
- RFID
- Fingerprint

Status and weighing history

- Ambient condition monitoring
- Digital balance auditor
- BWP guide
- Ambient light

Connectivity

- USB A + C, HDMI, Ethernet
- WiFi
- Hotspot (screen mirroring)

Applications

- | | |
|--------------------|-------------------------|
| ▪ Basic weighing | ▪ Pipette calibration |
| ▪ Parts counting | ▪ Differential weighing |
| ▪ Checkweighing | ▪ SQC |
| ▪ Dosing | ▪ Peak hold |
| ▪ Percent weighing | ▪ PGC |
| ▪ Density | ▪ Comparator |
| ▪ Animal weighing | ▪ Mass control |
| ▪ Formulations | ▪ Fraction |
| ▪ Statistics | |

Sartorius Cubis® II

Features and Selling Points

- More than 150 years market-proven expertise by an innovation leader in weighing
- Long lasting warranty with reliable service and global presence
- 7" Touch Display
- Automatic sliding doors with learning function
- Automatic Motorized Leveling with feet underneath the balance
- Ionizer
- isoCal
- Tool-free disassembly of draft shield and inner components of weighing compartment
- Titanium Weighing compartment
- Motion sensors for touch-free operation

User Log-In

- Password
- LDAP single sign on

Status and weighing history

- Status Center with most important information for simplified overview
- Ambient conditions can be monitored with a climate module

Connectivity

- USB A + B, Ethernet
- WiFi optional

Application Essentials

- | | |
|-----------------------|-------------------------------------|
| ▪ Averaging | ▪ Weighing |
| ▪ Multiplication | ▪ Totalization |
| ▪ Checkweighing | ▪ Interval print |
| ▪ Counting | ▪ Barcode print |
| ▪ Density of solids | ▪ Weighing with SBI output |
| ▪ Mixing | ▪ Additional applications available |
| ▪ Highest peak value | |
| ▪ Percentage weighing | |
| ▪ Statistics | |

Competitor Cross Guide: Ellipsis vs. Cubis® II

Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Minimum Weight (mg)*	Pan size (mm)	Chamber size DWH or DH (mm)
Ultramicro Balances									
UYA 2.5Y	0.0001	2.1	0.0015	0.0015	0.00015 0.00035	10 - 20	0.3	16	90 x 90
MCA2.7S-2S00-M	0.0001	2.1	0.0007	0.0005	0.00015 0.00018	7	0.3	20	80 x 67
UYA 2.5Y.F.A	0.0001	2.1	0.0015	0.0015	0.00015 0.00035	10 - 20	0.3	70 16	93.8 x 35
MCA2.7S-2S00-F	0.0001	2.1	0.0007	0.0005	0.00015 0.00018	7	0.3	50	109 x 33
Micro Balances									
MYA 2.5Y	0.001	2.1	0.003	0.003	0.00041 0.001	8	0.82	16	90 x 90
MYA 5.5Y	0.001	5.1	0.005	0.005	0.0006 0.001	8	1.2	26	90 x 90
MYA 6.5Y	0.001	6.1	0.005	0.005	0.0006 0.0016	8	1.2	26	90 x 90
MCA6.6S-2S00-M	0.001	6.1	0.003	0.003	0.0005 0.0006	5	0.82	30	80 x 67
MYA 5.5Y.F.A	0.001	5.1	0.005	0.005	0.0006 0.0016	8	1.2	70 16	93.8 x 35
MYA 5.5Y.F1	0.001	5.1	0.005	0.005	0.0006 0.0016	8	1.2	160 26	168 x 35
MCA6.6S-2S00-F	0.001	6.1	0.003	0.003	0.0005 0.0006	5	0.82	50	109 x 33
MYA 11.5Y	0.001	11	0.006	0.006	0.0009 0.0025	10	1.8	26	90 x 90
MCA10.6S-2S00-M	0.001	10.1	0.003	0.003	0.0005 0.0006	5	0.82	30	80 x 67
High Capacity Micro Balances									
XA 21.5Y.M.A	0.001	21	0.009	0.015	0.0013 0.0035	3.5	2.6	30	199 x 170 x 217
MCA36S-3S00-QP99-HWL	0.001	32	0.005	0.006	0.0007 0.0018	3.5	1.4	50 90	159 x 185 x 234
XA 52.5Y.M.A	0.005	52	0.02	0.02	0.0022 0.006	3.5	4.4	30	199 x 170 x 217
XA 53.5Y.M.A	0.001	53	0.02	0.02	0.0015 0.006	3.5	3	30	199 x 170 x 217
MCA66S-3S00-QP99-HWL	0.001	61	0.005	0.01	0.0007 0.0025	3.5	1.4	50 90	159 x 185 x 234
XA 6/21.5Y.M.A	0.001 0.002	6 21	0.009	0.015	0.0013 0.0035	3.5	2.6	30	199 x 170 x 217
MCA36P-3S00-QP99-HWL	0.001 0.01	10 32	0.006	0.008	0.0007 0.005	3.5 2.5	1.4	50 90	159 x 185 x 234
XA 21/52.5Y.M.A	0.001 0.005	21 52	0.02	0.02	0.0015 0.006	3.5	3	30	199 x 170 x 217
MCA66P-3S00-QP99-HWL	0.001 0.01	12 61	0.008	0.012	0.0007 0.006	3.5 2.5	1.4	50 90	159 x 185 x 234
Semi-Micro Balances									
XA 110.5Y.A	0.01	110	0.06	0.06	0.005 0.02	4	10	90 85	170 x 200 x 220
MCA125S-3S00-I-QP99	0.01	120	0.065	0.01	0.01 0.02	2	13	85 x 85	154 x 192 x 250
XA 210.5Y.A	0.01	210	0.1	0.1	0.005 0.025	4	10	90 85	170 x 200 x 220
MCA225S-3S00-I-QP99	0.01	220	0.065	0.01	0.01 0.02	2	13	85 x 85	154 x 192 x 250
XA 120 250.5Y.A	0.01 0.1	120 250	0.06	0.2	0.005 0.06	4	10	90 85	170 x 200 x 220
XA 82 220.5Y	0.01 0.1	82 220	0.06	0.2	0.005 0.06	4	10	90 85	170 x 200 x 220
MCA225P-3S00-D-QP99-HWL	0.01 0.1	120 220	0.03	0.06	0.007 0.02	1.5	13	90	185 x 159 x 234
Analytical Balances									
XA 220.5Y.A	0.1	220	0.2	0.2	0.05 0.08	1.3	100	100	170 x 200 x 220
MCA224S-2S00-I	0.1	220	0.13	0.12	0.05 0.05	1	100	85 x 85	154 x 192 x 250
XA 310.5Y.A	0.1	310	0.3	0.3	0.05 0.1	1.3	100	100	170 x 200 x 220
MCA324S-2S00-I	0.1	320	0.2	0.2	0.04 0.05	1	82	85 x 85	154 x 192 x 250
XA 520.5Y.A	0.1	520	0.5	0.4	0.07 0.18	1.3	140	100	170 x 200 x 220
MCA524S-2S00-I	0.1	520	0.2	0.2	0.04 0.05	1	82	85 x 85	154 x 192 x 250

Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Minimum Weight (mg)*	Pan size (mm)	Chamber size DWH or DH (mm)
Precision Balances									
PM 10.5Y	10	10,000	30	n/a	4 10	3		200 x 185	
MCA10202S-2S00-O	10	10,200	6	10	5 5	0.8		206 x 206	
PM 15.5Y	10	15,000	30	n/a	4 12	3		200 x 185	
MCA14202S-2S00-O	10	14,200	20	10	5 5	0.8		206 x 206	
High Capacity Balances									
PM 25.5Y	100	25,000	300	n/a	40 100	3		350 x 260	
PM 35.5Y	100	35,000	300	n/a	40 100	3		350 x 260	
MCA36201S-2S00-O	100	36,200	100	200	20 50	1.5		400 x 300	
PM 50.5Y	100	50,000	300	n/a	40 150	3		350 x 260	
MCA50201S-2S00-O	100	50,200	150	300	40 40	2		400 x 300	
PM 60.5Y	200	60,000	300	n/a	41 150	3		400 x 500	
MCA70201S-2S00-O	100	70,200	150	300	40 40	2		400 x 300	

* Typical values