

OPERATING SYSTEMS COMPATIBILITY CHART

COMPATIBILITY WITH MICROSOFT WINDOWS (JUNE 2023)

	Win 7 32 bit	Win 7 64 bit	Win 8 64 bit	Win 10 64 bit	Win 11 64 bit	remarks
MikroWin 2000	v.4.41	v.4.41	✗	✗	✗	
MikroWin 2010	✓	✓	✓*	✓*	✓*	* MikroWin 2010 v.5.21 or higher
ICE	✓	✓	✓	✓	✓*	* .Net Framework 2.0 required
ICE 2	✓	✓	✓	✓	✓	
LightCompass	✓	✓	✓	✓	✓	
Tubemaster	✓	✓	✓	✓	✗	Install as admin; run T_master.exe and Lcc.exe initially as admin
Winterm	✓	v.1.0.9	v.1.0.9	v.1.0.11	✗	For Berthold instruments, signed USB driver is needed
LBIS	v.3.20	v.3.20	v.4.0	v.4.0	v.4.0	
RadioStar v.4	✗	✗	✗	✗	✗	
RadioStar v.5	✓	✓	✓	✓	✓	
Indigo v.1.x	✓	✓	✓	✗	✗	Will not be supported anymore and can be replaced by Indigo v.2
Indigo v.2.x	✓	✓	✓	✓	✓	
Winglow	✓	✓	✗	✗	✗	

Version indicates the lowest compatible version

OPERATING SYSTEMS COMPATIBILITY CHART

COMPATIBILITY WITH MICROSOFT OFFICE (JUNE 2023)

	Office 2003	Office 2007	Office 2010	Office 2013	Office 2016	Office 365	remarks
MikroWin 2000							
MikroWin 2010							
ICE 2							
LightCompass							
Tubemaster		*	*	*	*	*	* Lumi.xls macro up to Excel 2003 only
Winterm							
LBIS							
RadioStar v.4	n.r.	n.r.	n.r.	n.r.	n.r.	n.r.	
RadioStar v.5	n.r.	n.r.	n.r.	n.r.	n.r.	n.r.	
Indigo v.1.x							
Indigo v.2.x							
Winglow							

n.r. = no relevance