

**Product** MicroBeta2  
**Model No's/part No** 2450/2460

产品中有害物质的名称及含量  
China EEP Hazardous Substances Information

部件名称 Component Name	有害物质 Hazardous Substances					
	铅	汞	镉	六价铬	多溴联苯	多溴二苯醚
	(Pb)	(Hg)	(Cd)	(Cr(VI))	(PBB)	(PBDE)
(电缆组件) Cable Assemblies	0	0	0	0	0	0
(显示器) Display	0	0	0	0	0	0
(电机组件) Electromechanical Assemblies	0	0	0	0	0	0
(紧固件) Fasteners	0	0	0	0	0	0
(机加工件) Machined Parts	0	0	0	0	0	0
(模具) Molded Parts	0	0	0	0	0	0
(电路板) PCBA's	0	0	0	0	0	0
(电源供应器) Power Supply	0	0	0	0	0	0
(钣金件) Sheetmetal	X	0	0	0	0	0
(管子) Tubing	0	0	0	0	0	0

本表格依据SJ/T11364的规定编制 This table is compiled according to SJ/T 11364 standard.

O: 表示该有害物质在该部件所有均质材料中的含量均在GB/T26572规定的限量要求以下。  
Indicates that the concentration of the hazardous substance in all homogeneous materials for the part is below the relevant threshold of the GB/T26572 standard.

X: 表示该有害物质至少在该部件的某一均质材料中的含量超出GB/T26572规定的限量要。  
Indicates that the concentration of the hazardous substance in at least one homogenous material of the part is above the relevant threshold of the GB/T 26572 standard.