

Product Wizard2 (IVD)
Model No's/part No 2470-0***/2480-0***/3470-0***

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	O	O	O	O	O	O
(显示器) Display	O	O	O	O	O	O
(电机组件) Electromechanical Assemblies	X	X	X	O	O	O
(紧固件) Fasteners	O	O	O	O	O	O
(机加工件) Machined Parts	O	O	O	O	O	O
(模具) Molded Parts	X	O	O	O	O	O
(电路板) PCBA's	O	O	O	O	O	O
(电源供应器) Power Supply	O	O	O	O	O	O
(钣金件) Sheetmetal	O	O	O	O	O	O
(管子) Tubing	O	O	O	O	O	O

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/T 26572 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.