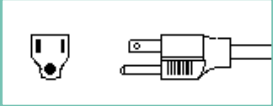
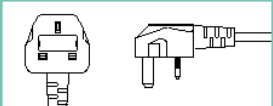
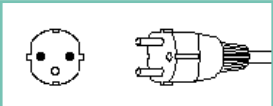
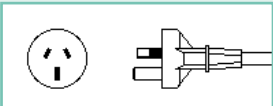
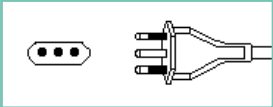
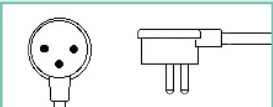



REGIONAL POWER CORDS

Description/Illustration	Catalogue Number
<p>EVOS® Power Cord, Type A (US)</p> <p>The Type A electrical plug is used in North America, Japan and parts of South America, Asia and the Middle East.</p> 	<p>AMEP4644</p>
<p>EVOS® Power Cord, Type G (UK)</p> <p>The Type G electrical plug is used in the UK, Ireland, Hong Kong, Singapore and other parts of Asia and Africa.</p> 	<p>AMEP4645</p>
<p>EVOS® Power Cord, Type C (EU)</p> <p>The Type C electrical plug is used throughout continental Europe, parts of the Middle East, and much of Africa, South America, Russia and the former Soviet republics.</p> 	<p>AMEP4646</p>
<p>EVOS® Power Cord, Type I (AU)</p> <p>The Type I electrical plug is mainly used in Australia and New Zealand, but also occurs in China and parts of Asia, South America and Africa.</p> 	<p>AMEP4647</p>
<p>EVOS® Power Cord, Type L (IT)</p> <p>The Type L plug is used in Italy, Chile, Uruguay and parts of North Africa.</p> 	<p>AMEP4648</p>
<p>EVOS® Power Cord, Type H (Israel)</p> <p>The Type H plug is used in Israel and Gaza.</p> 	<p>AMEP4649</p>
<p>EVOS® Power Cord, Type D (India/South Africa)</p> <p>The Type D plug is used in parts of Africa, Asia and South America.</p> 	<p>AMEP4708</p>