

**MicroTools™ Starter Kit 1**

MicroTools Kit 1 on 25 mm rods, preloaded into reusable magnetic bases

Cat. No.	Amount
T1SK-1	1 Kit

**Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** n/a**Content:**

- 20 Individual MicroTools on 25 mm rods, preloaded into reusable magnetic bases
- Magnetic Push-button Wand (for easily switching between tools)
- Serrated-end tweezers
- Holder and case
- Tool location placard

**nonsaleable to USA and Japan**