

## Incoterms 2000

### EXW

EX WORKS  
(...named place)

Ex works means that the seller delivers when he places the goods at the disposal of the buyer at the seller's premises or another named place ( i.e. works, factory, warehouse etc.) not cleared for export and not loaded on any collecting vehicle.

The term thus represents the minimum obligation for the seller, and the buyer has to bear all costs and risks involved in taking the goods from the seller's premises.