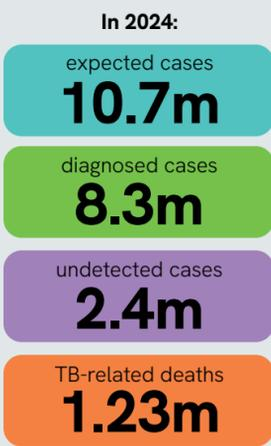
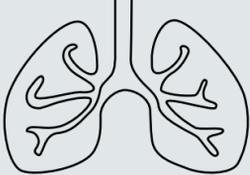
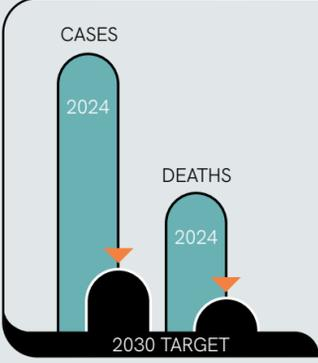


Join the fight against TB! Explore the challenging global TB landscape and the critical role of diagnostic testing in treatment, prevention and elimination.

## World TB headlines



TB is the **#1** infectious disease killer\*



TB remains a major global public health problem, with economic barriers restricting access to diagnosis and treatment. Although much progress is needed to meet the global target for 2030, incidence and deaths have reduced from 2023.

**Target:** The WHO's End TB strategy targets **80% drop** in new cases by 2030, **90% drop** in TB deaths

## A global health challenge

Nearly a **quarter of the global population** are estimated to be living with TB infection. While the majority of TB disease cases originate from a limited number of countries, **TB infection (LTBI) is widespread across the world.**

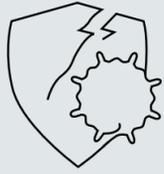


### At-risk populations and LTBI



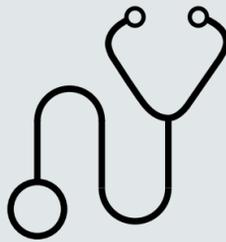
At-risk populations such as immunosuppressed, paediatric and HIV patients are most likely to progress from TB infection to TB disease - driving up fatalities and transmission. Highlighting this danger in 2024, **150,000 TB-related deaths occurred among people with HIV**, accounting for a significant proportion of TB mortality.

**Testing these populations is typically more difficult** due to low white blood cell counts.



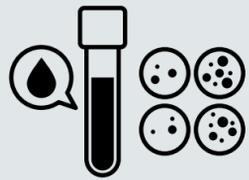
## Diagnosis is key

The TB infection reservoir is a silent driver of the epidemic - accurate and reliable diagnostic tests are key to breaking the cycle of transmission.

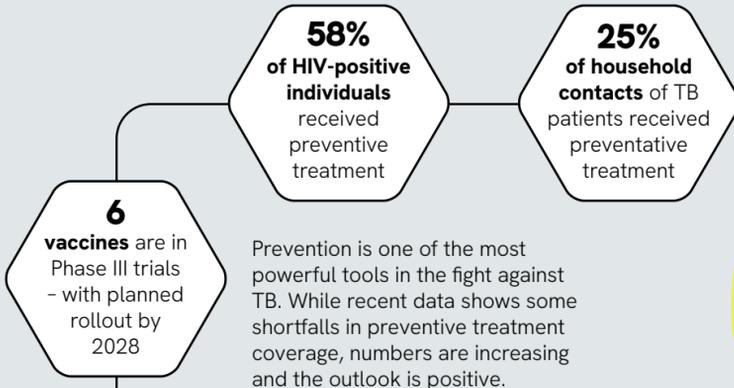


In 2024, the estimated prevalence of TB infection was about a quarter of the world's population.

**Increased use** of diagnostic tests is crucial to reducing LTBI, especially those that deliver **reliability, specificity and sensitivity in all patient groups** - including those most at risk.



## A path to prevention



Prevention is one of the most powerful tools in the fight against TB. While recent data shows some coverage, numbers are increasing and the outlook is positive.

**Target:** 90% preventative treatment coverage by 2027

## Taking the fight to TB

TB is once again the world's deadliest infectious disease, but **the path to elimination is clear:**



**Let's commit to amplifying awareness, driving innovation, and working together to achieve the WHO End TB Strategy's goals!**

