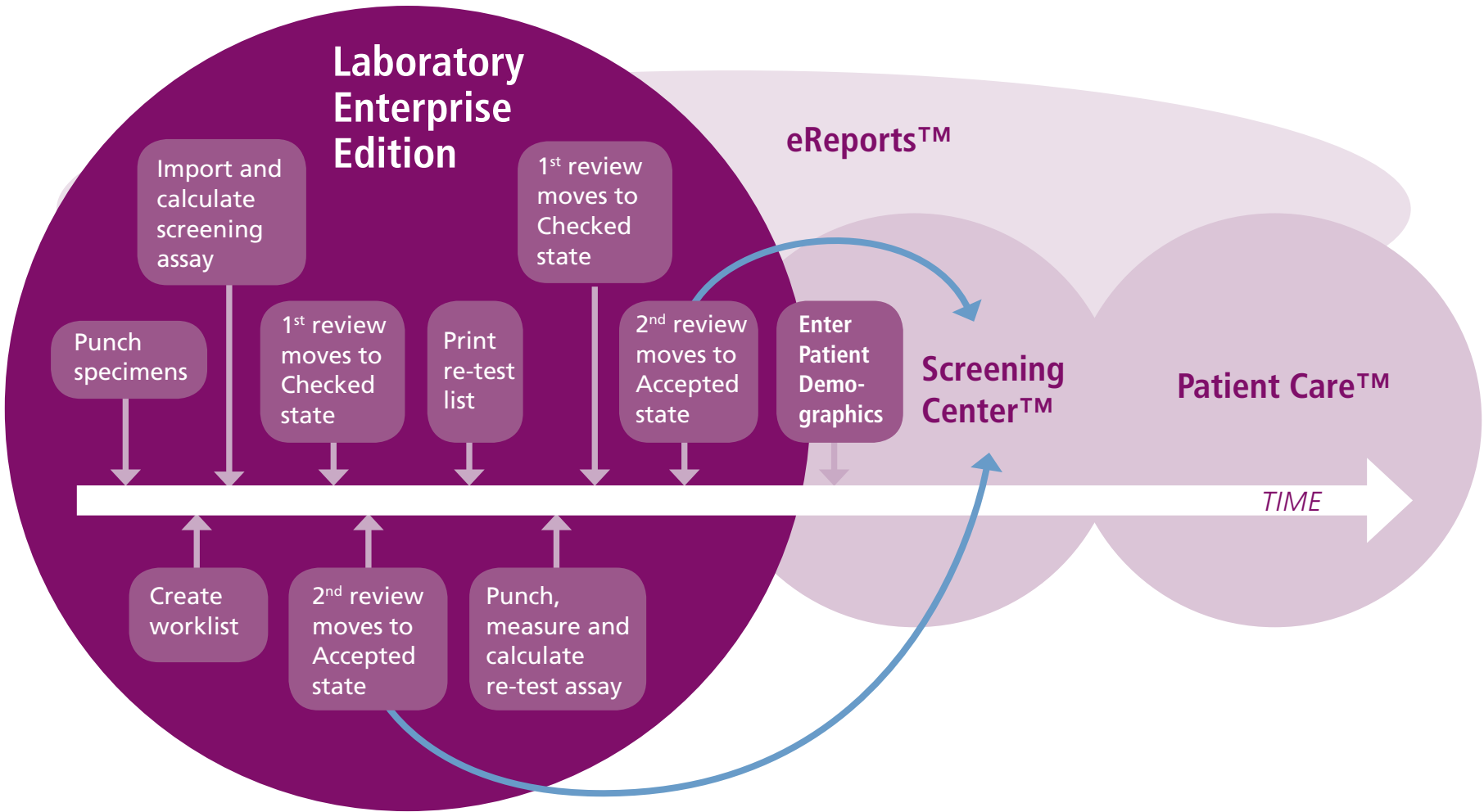
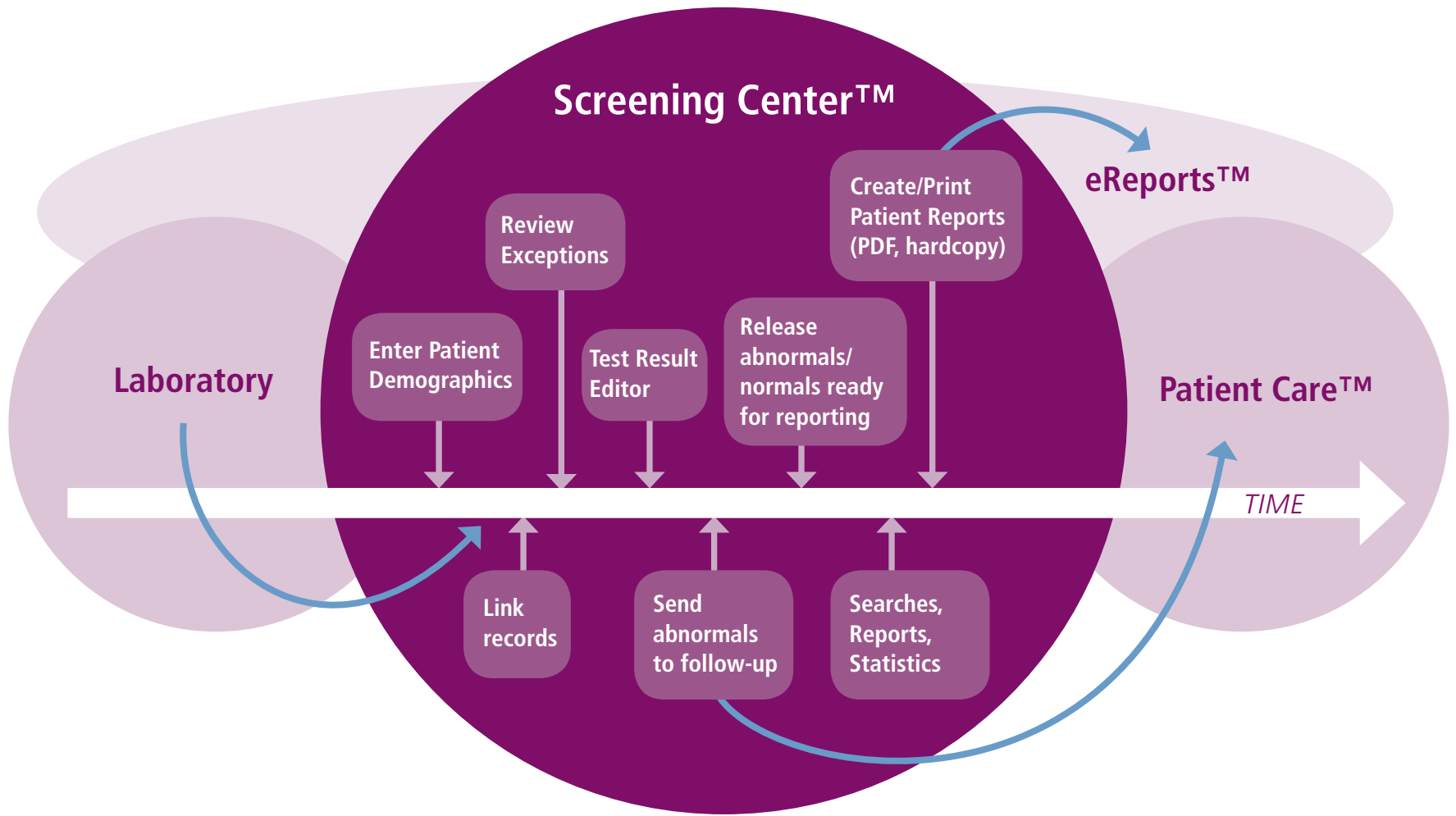


Specimen Gate® Data Flow



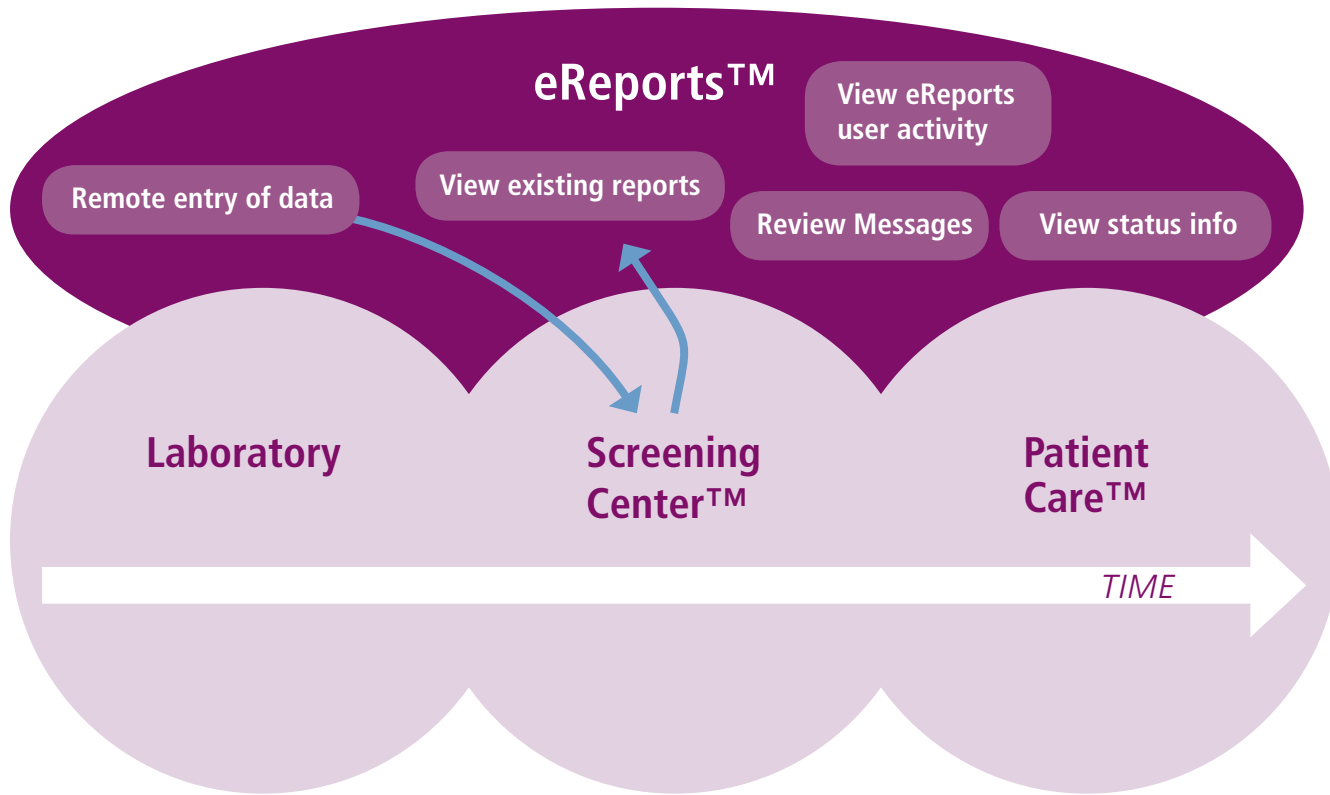
**Specimen Gate®
Data Flow**

Laboratory Enterprise Edition



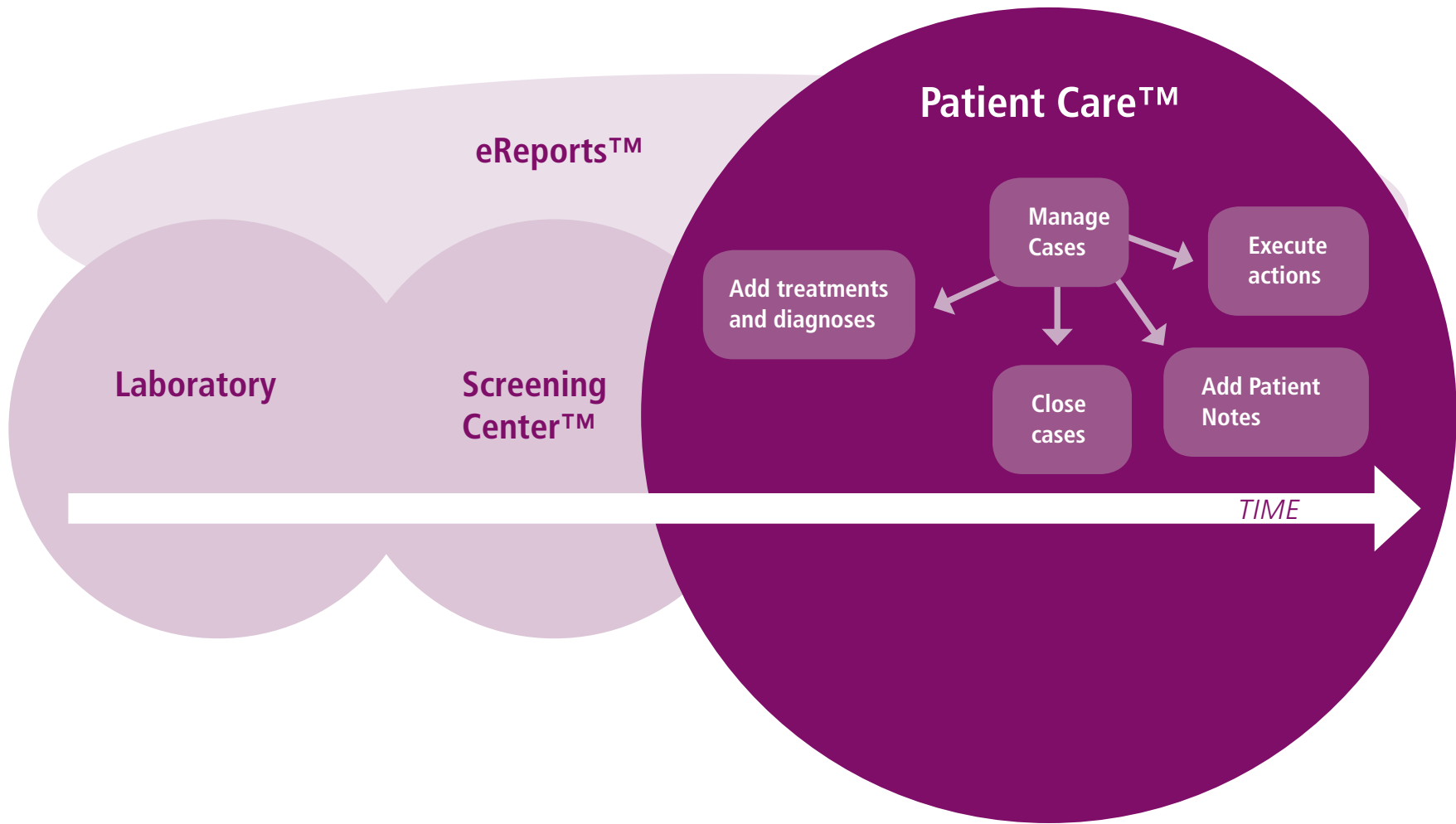
**Specimen Gate®
Data Flow**

Screening Center™



Specimen Gate® Data Flow

eReports™



Specimen Gate®
Data Flow

Patient Care™