

Real-time fluorescence assay for monitoring transglutaminase activity

Magdalena Adamczyk, Andreas Heil and Daniel Aeschlimann
Matrix Biology & Tissue Repair Research Unit, School of Dentistry, Cardiff University

- Isopeptidase assay for kinetic analysis of transglutaminase activity
- Optimized protocols allow for rapid implementation and standardization of measurements
- Assay is amenable to high throughput analysis of regulators/inhibitors of catalysis

Introduction

Transglutaminases (TGs) form a family of enzymes that catalyze various posttranslational protein modifications such as crosslinking, esterification and deamidation in a Ca^{2+} -dependent manner. Their main function is the formation of covalent N ϵ -[γ -glutamyl]lysine bonds within or between polypeptides to stabilize protein assemblies. The activity of these enzymes is crucial for tissue homeostasis and function in a number of organ systems, and the lack of or the excessive crosslinking activity have been linked to human disease processes. Here we perform kinetic measurements using recombinant TG2 and a fluorescent peptide model substrate in a format suitable for high-throughput analysis. This assay principle can be applied to kinetic studies on closely related enzymes including TG6 and can be optimised by modification of the backbone peptide sequence.

Assay Principle

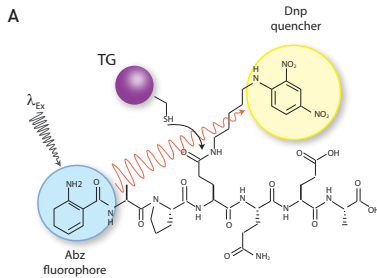


Fig. 1A: Structure of quenched substrate Abz-APE[γ -cad-Dnp]QEA whereby 2-aminobenzoyl (Abz) and 2,4-dinitrophenyl (Dnp) are fluorescent donor and quenching acceptor group, respectively.

The TG enzymatic reaction is a two-step process. The thioester intermediate of the enzyme formed with the substrate in the first step subsequently reacts with a nucleophile to regenerate active enzyme and release a 'crosslinked' polypeptide. The second step is reversible and, in the presence of an excess of crosslinked substrate, TG catalyzes isopeptide bond hydrolysis. We have exploited this latter activity for real-time monitoring of TG activity and characterize the effect of potential regulators on TG activity.

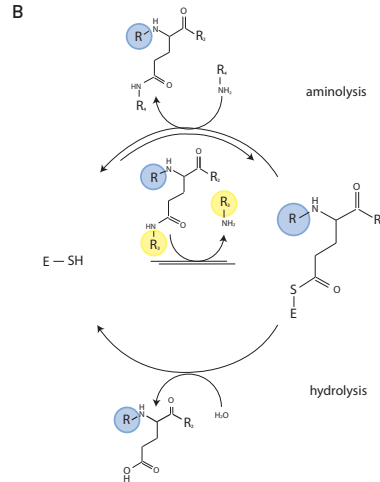


Fig. 1B: Reaction scheme for TG catalyzed isopeptidase reaction with substrate Abz-APE[γ -cad-Dnp]QEA in the presence of an excess of unlabeled primary amine [R'-NH $_2$, e.g. glycine methylester]. R: Abz-AP; R $_2$: QEA; R $_3$: [CH $_2$] $_n$ -NH-Dnp; E-SH [TG2 with active site thiol group].

Abz-APE[γ -cad-Dnp]QEA is a quenched fluorescent probe derived from a known glutamine donor substrate that mimics a crosslinked TG reaction product. In this peptide the fluorophore [2-aminobenzoyl (Abz)] is quenched by a 2,4-dinitrophenyl-cadaverine [cad-Dnp] substituent on the first Gln residue, essentially replacing the Lys side chain in N ϵ -[γ -glutamyl]lysine linked peptides (Fig. 1A). TG2-catalysed hydrolysis of the isopeptide bond releases the cad-Dnp moiety (Fig. 1B) and consequently generates an increase in light emission at $\lambda_{\text{max}}=418$ nm from the Abz group. The thioester enzyme intermediate formed is subsequently deacylated through either aminolysis or hydrolysis. Specificity of the reaction is guided by the amino acid residues that surround the reactive Gln residue.

Materials & Methods

- Microplate reader from BMG LABTECH
- Black optical bottom 96-well plates (Nunc)
- Abz-APE[γ -cad-Dnp]QEA TG2 substrate (Zedira). 50 mM stock in DMSO
- Transglutaminase 2, 1 mg/ml stock, from Zedira

Assay buffer

The assay buffer consists of 62.5 mM Tris/HCl, pH 7.4, 125 mM NaCl. Add glycine methylester (or alternative



amine donor substrate) and adjust pH immediately before use (at 37°C). Include DTT to prevent oxidative inactivation of TG.

Test protocol

Prime the microplate reader injectors with 20 mM CaCl₂ for enzyme activation (inj. 1) and H₂O or 20 mM MgCl₂ for control reaction (inj. 2). Pre-warm assay buffer and plate to 37°C and equilibrate instrument chamber at 37°C. Dilute substrate Abz-APE[γ-cad-Dnp]QEA (1:800) in assay buffer and add 80 µl of mixture into wells of the 96-well plate. Add desired amount of enzyme, e.g. 1 mg of TG2, and make up volume with H₂O to 90 µl. Transfer plate immediately into the reader and start program.

Reaction mixture (final concentrations)

The final assay volume is 100 µl and consists of 1-100 µg/ml TG2, 50 µM Abz-APE[γ-cad-Dnp]QEA, 10-55 mM glycine methylester or alternative nucleophile, as well as 1-5 mM DTT. After injection there is 2 mM CaCl₂ present in samples.

Instrument settings

Mode:	Fluorescence Intensity, plate mode
Filters:	Excitation: Ex320 Emission: 440-10
Optics:	top
No. of flashes:	20
Cycles:	90
Cycle time:	40 s (for 12 wells)
Injection cycle:	10
Injection volume:	10 µL
Shaking:	5 s after each cycle
Temperature:	37°C

TG2-mediated substrate conversion is linear for >30 min and initial reaction rates can be derived from linear regression of first 15-25 data points (Fig. 2A and B).

Conclusion

Reported here are optimized experimental conditions for determination of TG2 isopeptidase activity with the fluorescent model substrate Abz-APE[γ-cad-Dnp] QEA using the BMG LABTECH microplate readers to produce an assay that is rapid, direct and sensitive. Automated injection of Ca²⁺ for enzyme activation combined with the ability to continuously measure fluorescence intensity over a considerable time period with limited photo-bleaching facilitates the acquisition of kinetic data. Small sample size and plate format make the assay cost-effective and adaptable to high-throughput analysis.

Acknowledgements

This work was supported by grants from Coeliac UK and Arthritis Research UK (18461), and PhD studentships from Cardiff University/the President's Scholarship scheme to MA and AH.

Results & Discussion

After Ca²⁺ injection an increase in fluorescence can be observed dependent on the concentration of enzyme in sample (Fig. 2).

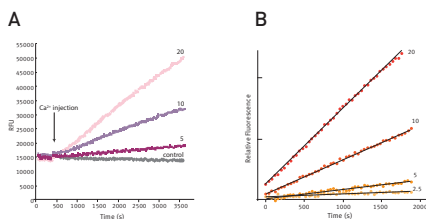


Fig. 2: Raw data [A] and processed data [B] of substrate Abz-APE[γ-cad-Dnp]QEA conversion at different concentrations of TG2 [2.5 – 20 µg/ml]. Control given represents 20 µg/ml TG2 without Ca²⁺ injection. Data processing involved normalization for well-specific fluorescence and subtraction of control to account for fluorescence bleaching.



PHERAstar® FSX

*The PHERAstar FSX is the newest PHERAstar reader.



CLARIOstar®



Omega Series