

Aβ1-42 binding and uptake by primary human astrocytes *in vitro*:

Effects of α1-Antichymotrypsin

H. M. Nielsen^{*1,2}, S. Janciauskiene², B. Holmqvist³, R. Veerhuis^{1,4}

*Researcher of the Alzheimer's Association; ICAAT Travel Fellowship 2006



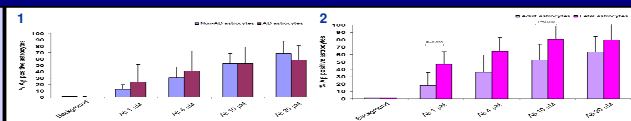
LUND UNIVERSITY
Faculty of Medicine

1) Departments of Clinical Chemistry, Pathology and the Alzheimer Center, VU Medical Center, Amsterdam, The Netherlands 2) Department of Clinical Sciences Malmö, Lund University, Malmö University

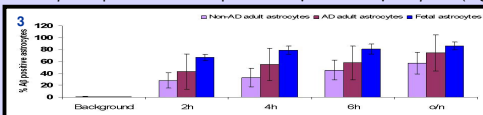
BACKGROUND

Imbalance between the production and clearance of the amyloid β-peptide (Aβ) is a key event in the Alzheimer's disease (AD) pathogenesis¹. Alpha1-antichymotrypsin (ACT) might influence the Aβ fibrillogenesis^{2,3}, biological effects^{4,5} and clearance⁶ leading to enhanced Aβ deposition in the brain⁷. Activated astrocytes are found surrounding amyloid plaques in AD brains⁸ but their role in AD pathogenesis is still poorly understood. These reactive cells over-express ACT⁹ and are able to release pro-inflammatory mediators¹⁰. Recent evidence also suggests that rodent astrocytes may internalize and degrade extracellular Aβ¹¹. If also human primary astrocytes are capable of degrading Aβ1-42, is still to be determined.

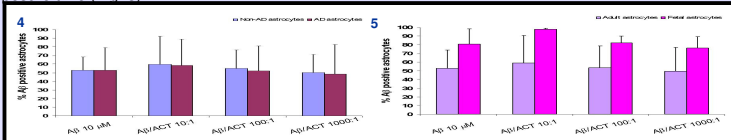
RESULTS



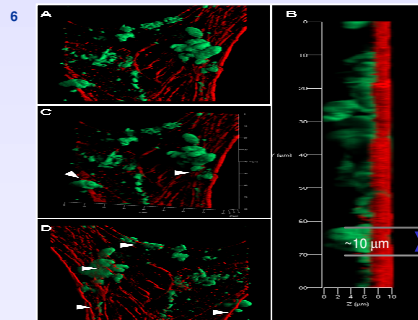
Non-AD (n=8) and AD-derived (n=6) adult astrocytes dose-dependently become Aβ1-42 positive in a similar manner upon o/n incubation (Fig. 1). Significantly more fetal astrocytes than adult astrocytes become Aβ1-42 positive when exposed to 1 μM and 10 μM Aβ1-42 (Fig. 2)



Adult non-AD (n=4), AD derived (n=3) and fetal astrocytes (n=4) become Aβ1-42 positive depending on exposure time (Fig. 3).



ACT in combination with Aβ1-42 has little or no effect on Aβ1-42 uptake by adult astrocytes (Fig. 4), whereas 17.34 % more fetal astrocytes become Aβ1-42 positive upon o/n treatment with 10 μM Aβ/ACT at a 10:1 molar ratio (Fig. 5)



Confocal Laser Scanning Microscopy optical sectioning (Z-scanning (10 μm stack) was performed on astrocytes incubated with FluoAβ (green) and labelled with TR conjugated phalloidin (F-actin, red) (Fig. 6).

Images show one Z-stack processed for 3D reconstruction. (A) View from above, FluoAβ aggregates of various size (green) are seen located on the cell surface, on top of the F-actin cytoskeleton.

(B) View (rotated 90°) of the astrocyte cell layer (10 μm) (red) where different size FluoAβ aggregates (green) are confirmed to preferentially be located on the cell surface (arrows).

(C) View diagonally from the side with FluoAβ aggregates seen intermingled with F-actin (arrows), indicating intracellular uptake.

(D) View from inside the cell (from below) showing various size FluoAβ aggregates (green) being intermingled with the F-actin (arrows).

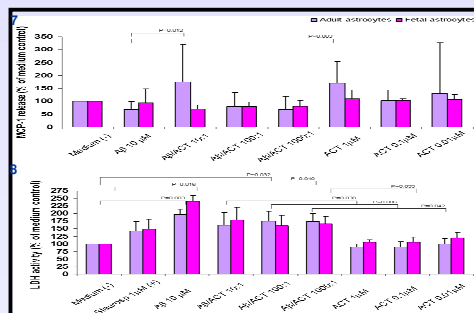
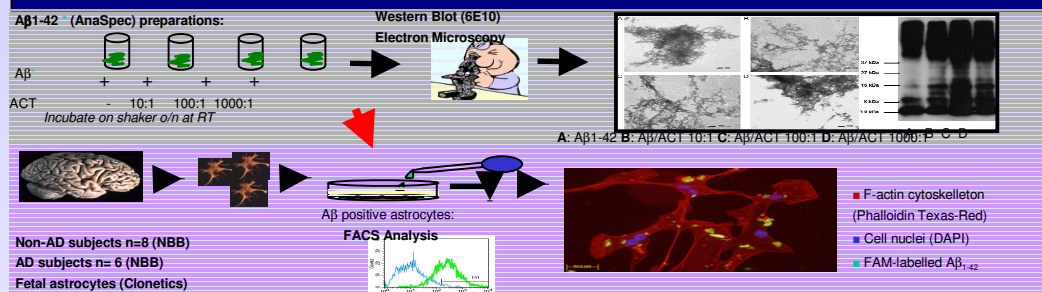
F-actin cytoskeleton (Phalloidin Texas-Red)

FAM-labelled Aβ₁₋₄₂

OBJECTIVE

- Can primary human fetal and adult astrocytes ingest Aβ1-42 *in vitro*?
- Do primary human fetal and adult astrocytes differ in *in vitro* Aβ1-42 uptake?
- Does α1-antichymotrypsin interfere with the Aβ uptake by primary human fetal and adult astrocytes?
- Is there a pro-inflammatory chemokine monocyte chemoattractant protein-1 (MCP-1) response to treatment with Aβ1-42 and Aβ/ACT combinations?

METHODS



The MCP-1 release from treated versus untreated cells was largely unaffected in both adult and fetal astrocytes (Fig. 7). However, in adult astrocytes the MCP-1 release was enhanced upon 10 μM Aβ/1 μM ACT (10:1) treatment versus 10 μM Aβ alone. This difference might be mediated by the effect of ACT itself.

Treatment with 10 μM Aβ o/n had a significant cytotoxic effect on both fetal and adult astrocytes (Fig. 8). In addition, the combinations of Aβ/ACT had a significantly more toxic effect on the astrocytes compared to the different concentrations of ACT alone.

CONCLUSIONS

- Primary human astrocytes are, during cytotoxic conditions, able to bind and take up Aβ1-42 *in vitro*, without a pro-inflammatory response
- Human adult astrocytes derived from non-AD and AD subjects become Aβ1-42 positive upon exposure, in a similar manner, whereas a greater percentage of human fetal astrocytes become Aβ1-42 positive upon 1 μM and 10 μM o/n treatment with Aβ1-42
- ACT has no or little effect on Aβ1-42 uptake by adult astrocytes whereas the uptake by fetal cells might be enhanced
- Enhanced MCP-1 release upon Aβ/ACT co-treatment versus Aβ alone in adult astrocytes, might be mediated by ACT itself